

**Important : Avant d'exercer votre choix, veuillez prendre connaissance des instructions situées au verso - Important : Before selecting please refer to instructions on reverse side**  
**Quelle que soit l'option choisie, noircir comme ceci ■ la ou les cases correspondantes, dater et signer au bas du formulaire - Whichever option is used, shade box(es) like this ■, date and sign at the bottom of the form**

☐ **JE DÉSIRE ASSISTER À CETTE ASSEMBLÉE** et demande une carte d'admission : dater et signer au bas du formulaire / **I WISH TO ATTEND THE SHAREHOLDER'S MEETING** and request an admission card : date and sign at the bottom of the form

## EUROFINS CEREP

S.A. à conseil d'administration  
au capital de 75 660 €  
Siège Social :  
Le Bois l'Evêque  
86600 Celle l'Evescault  
353 189 848 R.C.S. POITIERS

## ASSEMBLÉE GÉNÉRALE ORDINAIRE

Convoquée le 27 Mai 2021 à 11h00  
Au siège social, Le Bois l'Evêque 86600 Celle l'Evescault

### ORDINARY GENERAL MEETING

To be held on May 27th, 2021, at 11:00 a.m.  
At head office, Le Bois l'Evêque 86600 Celle l'Evescault

### CADRE RÉSERVÉ À LA SOCIÉTÉ - FOR COMPANY'S USE ONLY

Identifiant - Account

Vote simple  
Single vote

Vote double  
Double vote

Nominatif  
Registered

Porteur  
Bearer

Nombre d'actions  
Number of shares

Nombre de voix - Number of voting rights

|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------------------------|-----------------------------------------------------------------------------------------------------------|---|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|---|---|---|---|----|---|---|----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|-----------|--------------------------|------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|----------|--------------------------|--|--|--|--|--|--|--|--|--|--|------|--------------------------|----|----|----|----|----|----|----|----|----|----|---|---|----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|-----------|--------------------------|------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|----------|--------------------------|--|--|--|--|--|--|--|--|--|--|------|--------------------------|----|----|----|----|----|----|----|----|----|----|---|---|----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|-----------|--------------------------|------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|----------|--------------------------|--|--|--|--|--|--|--|--|--|--|------|--------------------------|----|----|----|----|----|----|----|----|----|----|---|---|----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|-----------|--------------------------|------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|----------|--------------------------|--|--|--|--|--|--|--|--|--|--|------|--------------------------|----|----|----|----|----|----|----|----|----|----|---|---|----------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|-----------|--------------------------|------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|--------------------------|----------|--------------------------|--|--|--|--|--|--|--|--|--|--|------|--------------------------|--|--|--|--|--|--|
| <input type="checkbox"/> <b>JE VOTE PAR CORRESPONDANCE / I VOTE BY POST</b><br>Cf. au verso (2) - See reverse (2)                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 |                          |                          |                          |                          |                          |                          |                          |                          |                          | Sur les projets de résolutions non agréés, je vote en noircissant la case correspondant à mon choix.<br>On the draft resolutions not approved, I cast my vote by shading the box of my choice. |                          | <input type="checkbox"/> <b>JE DONNE POUVOIR AU PRÉSIDENT DE L'ASSEMBLÉE GÉNÉRALE</b><br>Cf. au verso (3) |   | <input type="checkbox"/> <b>JE DONNE POUVOIR À :</b> Cf. au verso (4) pour me représenter à l'Assemblée<br><b>I HEREBY APPOINT:</b> See reverse (4) to represent me at the above mentioned Meeting<br>M. Mme ou Mlle, Raison Sociale / Mr, Mrs or Miss, Corporate Name |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Je vote <b>OUI</b> à tous les projets de résolutions présentés ou agréés par le Conseil d'Administration ou le Directoire ou la Gérance, à l'EXCEPTION de ceux que je signale en noircissant comme ceci ■ l'une des cases "Non" ou "Abstention". / I vote <b>YES</b> all the draft resolutions approved by the Board of Directors, EXCEPT those indicated by a shaded box, like this ■, for which I vote No or I abstain.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| <table border="1"> <tr> <td>1</td><td>2</td><td>3</td><td>4</td><td>5</td><td>6</td><td>7</td><td>8</td><td>9</td><td>10</td> <td>A</td><td>B</td> </tr> <tr> <td>Non / No</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Oui / Yes</td><td><input type="checkbox"/></td> </tr> <tr> <td>Abs.</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Non / No</td><td><input type="checkbox"/></td> </tr> <tr> <td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td> <td>Abs.</td><td><input type="checkbox"/></td> </tr> <tr> <td>11</td><td>12</td><td>13</td><td>14</td><td>15</td><td>16</td><td>17</td><td>18</td><td>19</td><td>20</td> <td>C</td><td>D</td> </tr> <tr> <td>Non / No</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Oui / Yes</td><td><input type="checkbox"/></td> </tr> <tr> <td>Abs.</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Non / No</td><td><input type="checkbox"/></td> </tr> <tr> <td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td> <td>Abs.</td><td><input type="checkbox"/></td> </tr> <tr> <td>21</td><td>22</td><td>23</td><td>24</td><td>25</td><td>26</td><td>27</td><td>28</td><td>29</td><td>30</td> <td>E</td><td>F</td> </tr> <tr> <td>Non / No</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Oui / Yes</td><td><input type="checkbox"/></td> </tr> <tr> <td>Abs.</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Non / No</td><td><input type="checkbox"/></td> </tr> <tr> <td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td> <td>Abs.</td><td><input type="checkbox"/></td> </tr> <tr> <td>31</td><td>32</td><td>33</td><td>34</td><td>35</td><td>36</td><td>37</td><td>38</td><td>39</td><td>40</td> <td>G</td><td>H</td> </tr> <tr> <td>Non / No</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Oui / Yes</td><td><input type="checkbox"/></td> </tr> <tr> <td>Abs.</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Non / No</td><td><input type="checkbox"/></td> </tr> <tr> <td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td> <td>Abs.</td><td><input type="checkbox"/></td> </tr> <tr> <td>41</td><td>42</td><td>43</td><td>44</td><td>45</td><td>46</td><td>47</td><td>48</td><td>49</td><td>50</td> <td>J</td><td>K</td> </tr> <tr> <td>Non / No</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Oui / Yes</td><td><input type="checkbox"/></td> </tr> <tr> <td>Abs.</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td> <td>Non / No</td><td><input type="checkbox"/></td> </tr> <tr> <td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td><td></td> <td>Abs.</td><td><input type="checkbox"/></td> </tr> </table> |                          |                          |                          |                          |                          |                          |                          |                          |                          | 1                                                                                                                                                                                              | 2                        | 3                                                                                                         | 4 | 5                                                                                                                                                                                                                                                                      | 6 | 7 | 8 | 9 | 10 | A | B | Non / No | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes | <input type="checkbox"/> | Abs. | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No | <input type="checkbox"/> |  |  |  |  |  |  |  |  |  |  | Abs. | <input type="checkbox"/> | 11 | 12 | 13 | 14 | 15 | 16 | 17 | 18 | 19 | 20 | C | D | Non / No | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes | <input type="checkbox"/> | Abs. | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No | <input type="checkbox"/> |  |  |  |  |  |  |  |  |  |  | Abs. | <input type="checkbox"/> | 21 | 22 | 23 | 24 | 25 | 26 | 27 | 28 | 29 | 30 | E | F | Non / No | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes | <input type="checkbox"/> | Abs. | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No | <input type="checkbox"/> |  |  |  |  |  |  |  |  |  |  | Abs. | <input type="checkbox"/> | 31 | 32 | 33 | 34 | 35 | 36 | 37 | 38 | 39 | 40 | G | H | Non / No | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes | <input type="checkbox"/> | Abs. | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No | <input type="checkbox"/> |  |  |  |  |  |  |  |  |  |  | Abs. | <input type="checkbox"/> | 41 | 42 | 43 | 44 | 45 | 46 | 47 | 48 | 49 | 50 | J | K | Non / No | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes | <input type="checkbox"/> | Abs. | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No | <input type="checkbox"/> |  |  |  |  |  |  |  |  |  |  | Abs. | <input type="checkbox"/> |  |  |  |  |  |  |
| 1                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 | 2                        | 3                        | 4                        | 5                        | 6                        | 7                        | 8                        | 9                        | 10                       | A                                                                                                                                                                                              | B                        |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Non / No                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes                                                                                                                                                                                      | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Abs.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                              | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No                                                                                                                                                                                       | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          | Abs.                                                                                                                                                                                           | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| 11                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                | 12                       | 13                       | 14                       | 15                       | 16                       | 17                       | 18                       | 19                       | 20                       | C                                                                                                                                                                                              | D                        |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Non / No                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes                                                                                                                                                                                      | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Abs.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                              | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No                                                                                                                                                                                       | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          | Abs.                                                                                                                                                                                           | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| 21                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                | 22                       | 23                       | 24                       | 25                       | 26                       | 27                       | 28                       | 29                       | 30                       | E                                                                                                                                                                                              | F                        |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Non / No                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes                                                                                                                                                                                      | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Abs.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                              | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No                                                                                                                                                                                       | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          | Abs.                                                                                                                                                                                           | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| 31                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                | 32                       | 33                       | 34                       | 35                       | 36                       | 37                       | 38                       | 39                       | 40                       | G                                                                                                                                                                                              | H                        |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Non / No                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes                                                                                                                                                                                      | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Abs.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                              | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No                                                                                                                                                                                       | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          | Abs.                                                                                                                                                                                           | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| 41                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                | 42                       | 43                       | 44                       | 45                       | 46                       | 47                       | 48                       | 49                       | 50                       | J                                                                                                                                                                                              | K                        |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Non / No                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Oui / Yes                                                                                                                                                                                      | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Abs.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                              | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | <input type="checkbox"/> | Non / No                                                                                                                                                                                       | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
|                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |                          |                          |                          |                          |                          |                          |                          |                          |                          | Abs.                                                                                                                                                                                           | <input type="checkbox"/> |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| Si des amendements ou des résolutions nouvelles étaient présentés en assemblée, je vote <b>NON</b> sauf si je signale un autre choix en noircissant la case correspondante :<br>In case amendments or new resolutions are proposed during the meeting, I vote <b>NO</b> unless I indicate another choice by shading the corresponding box:                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                        |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| - Je donne pouvoir au Président de l'assemblée générale. / I appoint the Chairman of the general meeting..... <input type="checkbox"/>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                            |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| - Je m'abstiens. / I abstain from voting ..... <input type="checkbox"/>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                           |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| - Je donne procuration [cf. au verso renvoi (4)] à M., Mme ou Mlle, Raison Sociale pour voter en mon nom ..... <input type="checkbox"/>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                           |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |
| I appoint [see reverse (4)] Mr, Mrs or Miss, Corporate Name to vote on my behalf..... <input type="checkbox"/>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                    |                          |                          |                          |                          |                          |                          |                          |                          |                          |                                                                                                                                                                                                |                          |                                                                                                           |   |                                                                                                                                                                                                                                                                        |   |   |   |   |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |    |    |    |    |    |    |    |    |    |    |   |   |          |                          |                          |                          |                          |                          |                          |                          |                          |                          |           |                          |      |                          |                          |                          |                          |                          |                          |                          |                          |                          |          |                          |  |  |  |  |  |  |  |  |  |  |      |                          |  |  |  |  |  |  |

Pour être pris en considération, tout formulaire doit parvenir au plus tard :

To be considered, this completed form must be returned no later than:

sur 1<sup>ère</sup> convocation / on 1st notification

sur 2<sup>ème</sup> convocation / on 2nd notification

à la banque / to the bank

22 mai 2021 / May 22nd, 2021

à la société / to the company

Date & Signature

« Si le formulaire est renvoyé daté et signé mais qu'aucun choix n'est coché (carte d'admission / vote par correspondance / pouvoir au président / pouvoir à mandataire), cela vaut automatiquement pouvoir au Président de l'assemblée générale »  
If the form is returned dated and signed but no choice is checked (admission card / postal vote / power of attorney to the President / power of attorney to a representative), this automatically applies as a proxy to the Chairman of the General Meeting

### 3) POUVOIR AU PRÉSIDENT DE L'ASSEMBLÉE GÉNÉRALE

Article L. 225-106 du Code de Commerce [extrait] :

\"Pour toute procuration d'un actionnaire sans indication de mandataire, le président de l'assemblée générale émet un vote favorable à l'adoption de projets de résolutions présentés ou agréés par le conseil d'administration ou le directoire, selon le cas, et un vote défavorable à l'adoption de tous les autres projets de résolution. Pour émettre tout autre vote, l'actionnaire doit faire choix d'un mandataire qui accepte de voter dans le sens indiqué par le mandant\".

à l'adoption de projets de résolutions présentés ou agréés par le conseil d'administration ou le directoire, selon le cas, et un vote défavorable à l'adoption de tous les autres projets de résolution. Pour émettre tout autre vote, l'actionnaire doit faire choix d'un mandataire qui accepte de voter dans le sens indiqué par le mandant\*.

## 23/12/2020