

Certificate TW17/00767.00 / 證書 TW17/00767.00

The management system of / 下述組織

SGS

Eurofins Panlabs Discovery Services Taiwan, Ltd 汎球生物科技股份有限公司

25, Wugong 6th Road, Wugu District, New Taipei City, 24891 Taiwan, R.O.C.
24891 新北市五股區五工六路 25 號

has been assessed and certified as meeting the requirements of
的管理系統已經過審核，並被證明符合下述要求

ISO 9001:2015

Quality management systems - Requirements

For the following activities / 所涵蓋的活動範圍

Provision of Nonclinical Drug Discovery and Development Solutions
非臨床醫藥之研發服務

This certificate is valid from 20 July 2026 until 20 July 2029 and remains valid subject to satisfactory surveillance audits.
此證書的有效期自 2026 年 07 月 20 日 至 2029 年 07 月 20 日 且其有效性應繫於持續符合的定期稽核

Issue 5. Certified since 20 July 2017

版次 5. 原始註冊日期 2017 年 07 月 20 日

Multiple certificates have been issued for this scope, the main certificate is numbered TW17/00767.00

在驗證範圍內出具了多張證書，主要證書編號為 TW17/00767.00

Certified activities performed by additional sites are listed on subsequent pages.

其它場區/合法實體之驗證活動於後續頁上列出

L. Moran

Authorised by / 簽署

Liz Moran

Business Manager

SGS United Kingdom Ltd

Rossmore Business Park, Ellesmere Port, Cheshire, CH65 3EN, UK

t +44 (0)151 350-6666 - www.sgs.com



This document is an authentic electronic certificate for Client' business purposes use only. Printed version of the electronic certificate are permitted and will be considered as a copy. This document is issued by the Company subject to SGS General Conditions of certification services available on [Terms and Conditions](#) | SGS. Attention is drawn to the limitation of liability, indemnification and jurisdictional clauses contained therein. This document is copyright protected and any unauthorized alteration, forgery or falsification of the content or appearance of this document is unlawful.
本文件為真實的電子版證書，僅供客戶用於其商業用途。此電子證書允許其列印版本並將被視為副本。本文件由本公司依據可取得之 [Terms and Conditions](#) | SGS。中之 SGS 驗證服務通用條款頒發，敬請注意其中亦已包含的責任範圍，賠償和司法管轄事項。本文件受版權保護，任何未經授權而對此文件的內容或外觀作出變更、偽造或篡改皆屬違法。



Eurofins Panlabs Discovery Services Taiwan, Ltd 汎球生物科技股份有限公司

ISO 9001:2015

Quality management systems - Requirements

Issue 5 / 版次 5

Sites / 場區

Pharmacology Discovery Services Taiwan, Ltd.

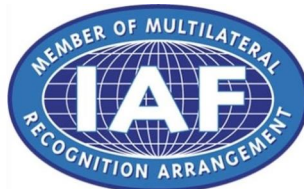
汎球藥理科技股份有限公司

25, Wugong 6th Road, Wugu District, New Taipei City, 24891 Taiwan, R.O.C.

24891 新北市五股區五工六路 25 號

Provision of Nonclinical Drug Discovery and Development Solutions

非臨床醫藥之研發服務



This document is an authentic electronic certificate for Client' business purposes use only. Printed version of the electronic certificate are permitted and will be considered as a copy. This document is issued by the Company subject to SGS General Conditions of certification services available on [Terms and Conditions](#) | SGS. Attention is drawn to the limitation of liability, indemnification and jurisdictional clauses contained therein. This document is copyright protected and any unauthorized alteration, forgery or falsification of the content or appearance of this document is unlawful.
本文件為真實的電子版證書，僅供客戶用於其商業用途。此電子證書允許其列印版本並將被視為副本。本文件由本公司依據可取得之 [Terms and Conditions](#) | SGS 中之 SGS 驗證服務通用條款頒發，敬請注意其中亦已包含的責任範圍，賠償和司法管轄事項。本文件受版權保護，任何未經授權而對此文件的內容或外觀作出變更、偽造或篡改皆屬違法。

