

Certificate SG23/00000060

The management system of

# Promega Pte Ltd



Business / Site Address: 77 Science Park Drive, #02-16, Cintech III Singapore Science Park 1, Singapore 118256

has been assessed and certified as meeting the requirements of  
**ISO 13485:2016**

For the following activities

1. Sales, storage, distribution of reagents and instruments systems used in In-vitro diagnostics.
2. Installation and servicing of instruments systems used in In-vitro diagnostics.

Technical Area:

Distribution Services, In Vitro Diagnostic Instruments and Software, Maintenance Services

This certificate is valid from 30 July 2024 until 11 October 2026 and remains valid subject to satisfactory surveillance audits.

Issue 2. Certified since 12 October 2023



Authorised by

Susan Law  
SEAP Regional Manager

SGS International Certification Services Singapore Pte Ltd  
30 Boon Lay Way, #03-01, Singapore 609957  
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.

