

Certificate SG23/00000036

The management system of

# Promega Pte Ltd

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**

**EN 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.

This certificate is valid from 15 October 2023 until 14 October 2026 and remains valid subject to satisfactory surveillance audits.

Issue 2. Certified since 26 July 2023

Organization certified since 15 December 2008 and first certified by SGS on 26 July 2023.

Authorised by  
Jonathan Hall  
Global Head - Certification Services

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.

