

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.

This certificate is valid from 12 October 2023 until 14 October 2026 and remains valid subject to satisfactory surveillance audits.

Issue 1. 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.

