

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Promega Pte Ltd
77 Science Park Drive
#02-16
Cintech III Singapore Science Park 1
Singapore 118256

Holds Certificate No:

MD 541260

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 for the following scope:

The provision of sales & distribution of products, service and information applicable to the delivery of reagents, reagents support systems, instrumentation and medical devices for the biotechnology, DNA identification and in vitro diagnostic markets

For and on behalf of BSI:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2008-12-15

Latest Revision Date: 2020-09-18

Effective Date: 2020-10-15

Expiry Date: 2023-10-14

Page: 1 of 1



...making excellence a habit.™

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsi-global.com/ClientDirectory or telephone +(65)62700777.

Further clarifications regarding the scope of this certificate and the applicability of ISO 13485:2016 requirements may be obtained by consulting the organization.

This certificate is valid only if provided original copies are in complete set.