

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Promega Biosystems Inc.
1017,A Building Gangseo-Hangang-Xi Tower
401, Yangcheon-ro
Gangseo-gu
Seoul
07528
Republic of Korea

Holds Certificate Number: MD 558018

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

The design and manufacture of Nucleic Acid and Protein auto extraction devices.

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Global Regulatory & Quality

Original Registration Date: 2010-01-20

Latest Revision Date: 2024-11-29

Effective Date: 2025-01-08

Expiry Date: 2028-01-07

Page: 1 of 1



...making excellence a habit.™