

# 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:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2010-01-20

Latest Revision Date: 2021-12-30

Effective Date: 2022-01-08

Expiry Date: 2025-01-07

Page: 1 of 1



...making excellence a habit.™