

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 No:

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:



Stewart Brain, Head of Compliance & Risk - Medical Devices

Original Registration Date: 2010-01-20

Latest Revision Date: 2018-12-24

Effective Date: 2019-01-08

Expiry Date: 2022-01-07

Page: 1 of 1



003

...making excellence a habit.™