

# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

**Promega Corporation**  
2800 Woods Hollow Road  
Madison  
Wisconsin  
53711-5399  
USA

Holds Certificate Number:

FM 505517

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

The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2006-08-24

Latest Revision Date: 2023-02-08

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 1 of 3



...making excellence a habit.™

Certificate No: FM 505517

| Location   | Registered Activities   |
|--|---|
| Promega Corporation<br>RDC<br>2800 Woods Hollow Road<br>Madison<br>Wisconsin<br>53711-5399<br>USA        | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>Feynman Center<br>2780 Woods Hollow Drive<br>Madison<br>Wisconsin<br>53711<br>USA | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>Kepler Center<br>5455 Nobel Drive<br>Madison<br>Wisconsin<br>53711<br>USA         | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>BTC<br>5445 East Cheryl Parkway<br>Madison<br>Wisconsin<br>53711<br>USA           | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>RFC<br>2617 Progress Rd<br>Madison<br>Wisconsin<br>53716<br>USA                   | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |

Original Registration Date: 2006-08-24

Latest Revision Date: 2023-02-08

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 2 of 3

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.bsigroup.com/ClientDirectory](http://www.bsigroup.com/ClientDirectory)

Issuing Body: BSI Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.

Contact Office: 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA.

Certificate No: FM 505517

| Location   | Registered Activities   |
|--|---|
| Promega Corporation<br>Kornberg Center<br>5430 East Cheryl Parkway<br>Madison<br>Wisconsin<br>53711<br>USA           | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>Instrumentation Building<br>5500 & 5510 Nobel Dr.<br>Madison<br>Wisconsin<br>53711<br>USA     | The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets. |
| Promega Corporation<br>Chapelle Manufacturing Center<br>3075 Sub-Zero Pkwy<br>Fitchburg<br>Wisconsin<br>53719<br>USA | Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.   |
| Promega Corporation<br>Faraday<br>2800 Fish Hatchery Road<br>Madison<br>Wisconsin<br>53711<br>USA                    | Manufacturing.  |

Original Registration Date: 2006-08-24

Latest Revision Date: 2023-02-08

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 3 of 3

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.bsigroup.com/ClientDirectory](http://www.bsigroup.com/ClientDirectory)

Issuing Body: BSI Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.

Contact Office: 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA.