

# 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

Facility ID Number: F005139

Holds Certificate No:

**MDSAP 732547**

The company listed on this certificate has been audited to and found to conform with ISO 13485:2016 including the following country specific requirements:

**Canada:** Medical Devices Regulations - Part 1 - SOR 98/282

**USA:** 21 CFR 820, 21 CFR 803, 21 CFR 806, 21 CFR 807 - Subparts A to D

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

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Global Regulatory & Quality

Original Registration Date: 2021-08-11

Effective Date: 2025-10-22

Expiry Date: 2027-08-08



BSI Group America Inc. is an MDSAP recognised auditing organization

Page: 1 of 4

...making excellence a habit.™

Certificate No: **MDSAP 732547**

Location	Registered Activities
Promega Corporation 2800 Woods Hollow Road Madison Wisconsin 53711-5399 USA Facility ID Number: F005139	Fermentation and Manufacturing.
Promega Corporation 2780 Woods Hollow Drive Madison Wisconsin 53711 USA Facility ID Number: F005139	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 5455 Nobel Drive Madison Wisconsin 53711 USA Facility ID Number: F005139	Receiving, Manufacturing, Packaging, Shipping, Warehouse.
Promega Corporation 5445 East Cheryl Parkway Madison Wisconsin 53711 USA Facility ID Number: F005139	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 2617 Progress Rd Madison Wisconsin 53716 USA Facility ID Number: F005139	Receiving, Manufacturing, Warehouse.

Original Registration Date: 2021-08-11

Effective Date: 2025-10-22

Expiry Date: 2027-08-08

Page: 2 of 4

This certificate remains the property of BSI and shall be returned immediately upon request.  
An electronic certificate can be authenticated [online](https://www.bsigroup.com/ClientDirectory). Printed copies can be validated at [www.bsigroup.com/ClientDirectory](https://www.bsigroup.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.

Americas Headquarters: BSI Group America Inc., 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA  
A Member of the BSI Group of Companies.

Certificate No: **MDSAP 732547**

Location	Registered Activities
Promega Corporation 5500 East Cheryl Parkway Madison Wisconsin 53711 USA Facility ID Number: F005139	Administrative Offices.
Promega Corporation 2798 South Fish Hatchery Road Madison Wisconsin 53711 USA Facility ID Number: F005139	Engineering & Operations Support.
Promega Corporation 5430 East Cheryl Parkway Madison Wisconsin 53711 USA Facility ID Number: F005139	Research and Development, Manufacturing.
Promega Corporation 5511 Bjorksten Place Madison Wisconsin 53711 USA Facility ID Number: F005139	Engineering & Operations Support.
Promega Corporation 5500/5510 Nobel Drive Madison Wisconsin 53711 USA Facility ID Number: F005139	Research and Development, Receiving, Manufacturing, Packaging, Warehouse, Instrument Servicing.

Original Registration Date: 2021-08-11

Effective Date: 2025-10-22

Expiry Date: 2027-08-08

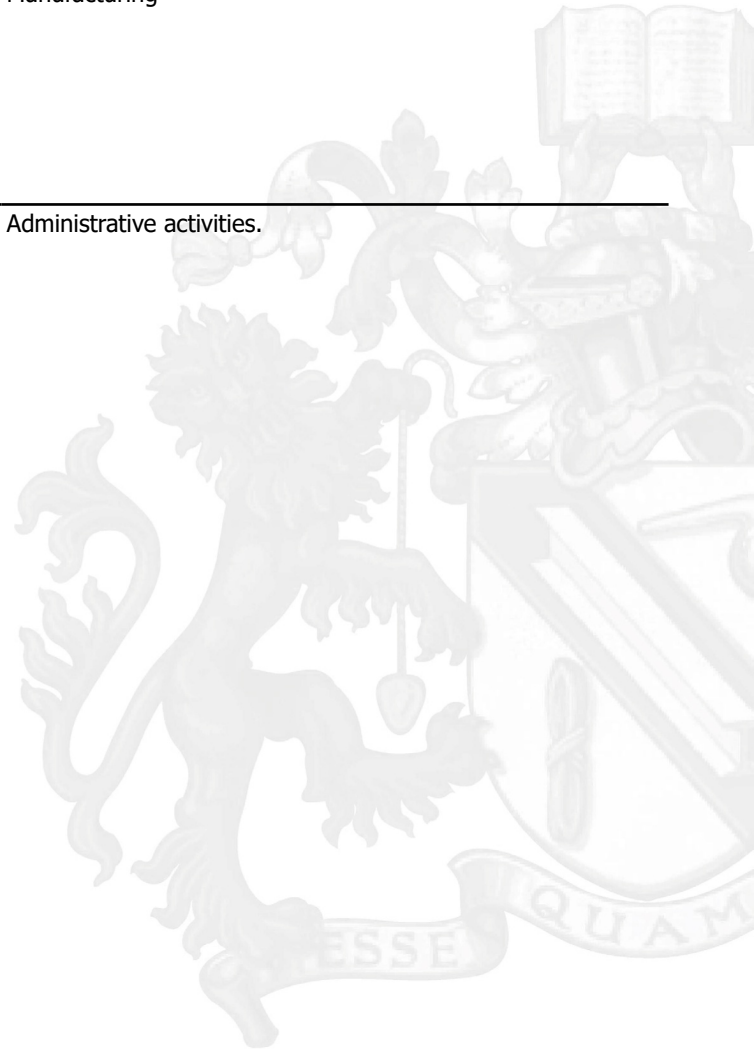
Page: 3 of 4

This certificate remains the property of BSI and shall be returned immediately upon request.  
An electronic certificate can be authenticated [online](https://www.bsigroup.com/ClientDirectory). Printed copies can be validated at [www.bsigroup.com/ClientDirectory](https://www.bsigroup.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.

Americas Headquarters: BSI Group America Inc., 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA  
A Member of the BSI Group of Companies.

Certificate No: **MDSAP 732547**

Location	Registered Activities
Promega Corporation 3075 Sub-Zero Pkwy Fitchburg Wisconsin 53719 USA Facility ID Number: F005139	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 2800 Fish Hatchery Road Madison Wisconsin 53711 USA Facility ID Number: F005139	Manufacturing
Promega Corporation 5520 Research Park Drive Madison Wisconsin 53711 USA Facility ID Number: F005139	Administrative activities.



Original Registration Date: 2021-08-11

Effective Date: 2025-10-22

Expiry Date: 2027-08-08

Page: 4 of 4

This certificate remains the property of BSI and shall be returned immediately upon request.  
An electronic certificate can be authenticated [online](https://www.bsigroup.com/ClientDirectory). Printed copies can be validated at [www.bsigroup.com/ClientDirectory](https://www.bsigroup.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.

Americas Headquarters: BSI Group America Inc., 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA  
A Member of the BSI Group of Companies.