

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

Statement of Conformity: The company listed on this certificate has been audited to and found to conform with the following criteria: ISO 13485:2016 and 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:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2021-08-11

Effective Date: 2021-08-11

Expiry Date: 2024-08-08



BSI Group America Inc. is an MDSAP authorized auditing organization

Page: 1 of 3

...making excellence a habit.™

Certificate No: **MDSAP 732547**

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

Original Registration Date: 2021-08-11

Effective Date: 2021-08-11

Expiry Date: 2024-08-08

Page: 2 of 3

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
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 Agora 5500 East Cheryl Parkway Madison Wisconsin 53711 USA Facility ID Number: F005139	Administrative Offices.
Promega Corporation Lamarr Center 2798 South Fish Hatchery Road Madison Wisconsin 53711 USA Facility ID Number: F005139	Engineering & Operations Support.
Promega Corporation Kornberg Center 5430 East Cheryl Parkway Madison Wisconsin 53711 USA Facility ID Number: F005139	Research and Development.
Promega Corporation daVinci 5511 Bjorksten Place Madison Wisconsin 53711 USA Facility ID Number: F005139	Engineering & Operations Support.
Promega Corporation Instrumentation Building 5500 Nobel Drive Madison Wisconsin 53711 USA Facility ID Number: F005139	Manufacturing of instruments, technical services, and the installation and servicing of all instruments.

Original Registration Date: 2021-08-11

Effective Date: 2021-08-11

Expiry Date: 2024-08-08

Page: 3 of 3

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
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.