

# 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 No:

**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:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2006-08-24

Latest Revision Date: 2021-08-11

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 RDC 2800 Woods Hollow Road Madison Wisconsin 53711-5399 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.
Promega Corporation Feynman Center 2780 Woods Hollow Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.
Promega Corporation Kepler Center 5455 Nobel Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.
Promega Corporation BTC 5445 East Cheryl Parkway Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.
Promega Corporation RFC 2617 Progress Rd Madison Wisconsin 53716 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: 2021-08-11

Effective Date: 2021-08-09

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](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: **FM 505517**

Location

Promega Corporation  
Kornberg Center  
5430 East Cheryl Parkway  
Madison  
Wisconsin  
53711  
USA

Registered Activities

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

Promega Corporation  
Instrumentation Building  
5500 Nobel Drive  
Madison  
Wisconsin  
53711  
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: 2021-08-11

Effective Date: 2021-08-09

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](#). Printed copies can be validated at [www.bsigroup.com/ClientDirectory](http://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.