



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:


Chief Operating Officer Assurance - Americas

Original Registration Date: 2006-08-24

Latest Revision Date: 2018-06-13

Effective Date: 2018-08-24

Expiry Date: 2021-08-23

Page: 1 of 2



...making excellence a habit.™

Certificate No: **FM 505517**

Location	Registered Activities
Promega Corporation 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 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 8137 Forsythia St. Middleton Wisconsin 53562 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation for the medical device and in-vitro diagnostics markets.
Promega Corporation 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 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 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

Effective Date: 2018-08-24

Latest Revision Date: 2018-06-13

Expiry Date: 2021-08-23

Page: 2 of 2

This certificate remains the property of BSI and shall be returned immediately upon request.
An electronic certificate can be authenticated [online](http://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.