



Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

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

Holds Certificate No:

FM 591295

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

The design, manufacture and servicing of reagents, reagent support systems and instrumentation.

For and on behalf of BSI:


Chief Operating Officer Assurance - Americas

Original Registration Date: 2012-09-13

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 591295**

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.
Promega Corporation 2780 Woods Hollow Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation 8137 Forsythia St. Middleton Wisconsin 53562 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation 5455 Nobel Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation 5445 East Cheryl Parkway Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation 2617 Progress Rd Madison Wisconsin 53716 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.

Original Registration Date: 2012-09-13

Effective Date: 2018-08-24

Latest Revision Date: 2018-06-13

Expiry Date: 2021-08-23

Page: 2 of 2