

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:

Carlos Pitanga, Chief Operating Officer Assurance – Americas

Original Registration Date: 2012-09-13

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 591295**

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.
Promega Corporation Feynman Center 2780 Woods Hollow Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation Kepler Center 5455 Nobel Drive Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation BTC 5445 East Cheryl Parkway Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation RFC 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

Latest Revision Date: 2021-08-11

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 2 of 3

Certificate No: **FM 591295**

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.

Promega Corporation
Instrumentation Building
5500 Nobel Drive
Madison
Wisconsin
53711
USA

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



Original Registration Date: 2012-09-13

Latest Revision Date: 2021-08-11

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 3 of 3