

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 Number:

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:

Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2006-08-24

Latest Revision Date: 2023-08-31

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: 2023-08-31

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 2 of 3

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsigroup.com/ClientDirectory

Issuing Body: BSI Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.

Contact Office: 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA.

Certificate No: FM 505517

Location	Registered Activities
Promega Corporation Kornberg Center 5430 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 Alan Turing Center (ATC) 5500/5510 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 Chappelle Manufacturing Center 3075 Sub-Zero Pkwy Fitchburg Wisconsin 53719 USA	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation Faraday 2800 Fish Hatchery Road Madison Wisconsin 53711 USA	Manufacturing.

Original Registration Date: 2006-08-24

Latest Revision Date: 2023-08-31

Effective Date: 2021-08-09

Expiry Date: 2024-08-08

Page: 3 of 3

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsigroup.com/ClientDirectory

Issuing Body: BSI Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.

Contact Office: 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA.