

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 diagnostic markets.

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Global Regulatory & Quality

Original Registration Date: 2006-08-24

Latest Revision Date: 2024-08-02

Effective Date: 2024-08-09

Expiry Date: 2027-08-08

Page: 1 of 3



...making excellence a habit.™

Certificate No: FM 505517

Location	Registered Activities
Promega Corporation 2800 Woods Hollow Road Madison Wisconsin 53711-5399 USA	Fermentation and Manufacturing.
Promega Corporation 2780 Woods Hollow Drive Madison Wisconsin 53711 USA	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 5455 Nobel Drive Madison Wisconsin 53711 USA	Receiving, Manufacturing, Packaging, Shipping, Warehouse.
Promega Corporation 5445 East Cheryl Parkway Madison Wisconsin 53711 USA	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 2617 Progress Rd Madison Wisconsin 53716 USA	Receiving, Manufacturing, Warehouse.
Promega Corporation 5500 East Cheryl Parkway Madison Wisconsin 53711 USA	Administrative Offices.

Original Registration Date: 2006-08-24

Latest Revision Date: 2024-08-02

Effective Date: 2024-08-09

Expiry Date: 2027-08-08

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 2798 South Fish Hatchery Road Madison Wisconsin 53711 USA	Engineering & Operations Support.
Promega Corporation 5430 East Cheryl Parkway Madison Wisconsin 53711 USA	Research and Development, Manufacturing.
Promega Corporation 5511 Bjorksten Place Madison Wisconsin 53711 USA	Engineering & Operations Support.
Promega Corporation 5500/5510 Nobel Drive Madison Wisconsin 53711 USA	Research and Development, Receiving, Manufacturing, Packaging, Warehouse, Instrument Servicing.
Promega Corporation 3075 Sub-Zero Pkwy Fitchburg Wisconsin 53719 USA	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.
Promega Corporation 2800 Fish Hatchery Road Madison Wisconsin 53711 USA	Manufacturing.

Original Registration Date: 2006-08-24

Latest Revision Date: 2024-08-02

Effective Date: 2024-08-09

Expiry Date: 2027-08-08

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.