

Certificate of Registration

ENVIRONMENTAL MANAGEMENT SYSTEM - ISO 14001:2015

This is to certify that:

Promega Corporation
Arnold Center
2800 Woods Hollow Road
Madison
Wisconsin
53711
USA

Holds Certificate No:

EMS 805637

and operates an Environmental Management System which complies with the requirements of ISO 14001:2015 for the following scope:

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

For and on behalf of BSI:

Jessica Patel, Senior Vice President, Assurance Americas

Original Registration Date: 2025-04-16

Effective Date: 2025-04-16

Latest Revision Date: 2025-04-16

Expiry Date: 2028-04-15

Page: 1 of 4



...making excellence a habit.™

Certificate No: **EMS 805637**

Location	Registered Activities
Promega Corporation Arnold Center 2800 Woods Hollow Road Madison Wisconsin 53711 USA	Administrative Offices, Wellness Center.
Promega Corporation Agora 5500 East Cheryl Parkway Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
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.
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 Chappelle Manufacturing Center 3075 Sub-Zero Pkwy Fitchburg Wisconsin 53719 USA	Receiving, Manufacturing, Dispensing, Packaging, Shipping, Warehouse.

Original Registration Date: 2025-04-16

Latest Revision Date: 2025-04-16

Effective Date: 2025-04-16

Expiry Date: 2028-04-15

Page: 2 of 4

Certificate No: **EMS 805637**

Location	Registered Activities
Promega Corporation 2800 Fish Hatchery Road Madison Wisconsin 53711 USA	Manufacturing.
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 Hangar 3792 Corbin Court Madison Wisconsin 53704 USA	Airplane Hangar.
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 Kornberg Center 5430 East Cheryl Parkway Madison Wisconsin 53711 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.

Original Registration Date: 2025-04-16

Latest Revision Date: 2025-04-16

Effective Date: 2025-04-16

Expiry Date: 2028-04-15

Page: 3 of 4

Certificate No: **EMS 805637**

Location	Registered Activities
Promega Corporation RFC 2617 Progress Rd Madison Wisconsin 53716 USA	The design, manufacture and servicing of reagents, reagent support systems and instrumentation.
Promega Corporation daVinci Center 5511 Bjorksten Place Madison Wisconsin 53711 USA	Operations Support, Engineering, Office.
Promega Corporation Lamarr Center 2798 South Fish Hatchery Road Madison Wisconsin 53711 USA	Operations Support, Engineering, Office.

Original Registration Date: 2025-04-16

Latest Revision Date: 2025-04-16

Effective Date: 2025-04-16

Expiry Date: 2028-04-15

Page: 4 of 4