

CERTIFICATE OF REGISTRATION

This is to certify that the management system of:

Promega Corporation

Main Site: 277 Granada Drive, San Luis Obispo,
California 93401, United States

has been registered by Intertek as conforming to the requirements of:

ISO 9001:2015

The management system is applicable to:

Design and Manufacture of Specialty Chemicals.

Certificate Number:

0154719

Initial Certification Date:

05 April 2005

Last Certificate Expiry Date:

07 August 2023

Date of Last Recertification Audit:

28 June 2023

Certification Cycle Start Date:

11 August 2023

Issuing Date:

11 August 2023

Valid Until:

07 August 2026



intertek

MANAGEMENT SYSTEMS
CERTIFICATION BODY

Calin Moldovean

President, Business Assurance

Intertek Testing Services NA, Inc. 900 Chelmsford
Street, Lowell MA 01851, USA

