

# 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

**Date of Certification Decision:**

14 June 2026

**Issuing Date:**

14 June 2026

**Valid Until:**

07 August 2029



intertek



A handwritten signature in black ink that reads "Mark Thomas".

**Mark Thomas**

Executive Vice President

Assurance, Sustainability, AgriWorld &  
Food

Intertek Testing Services NA,  
Inc. dba Intertek,  
4700 Broadmoor SE Suite 200 Kentwood,  
Michigan 49512, United States

