



# CERTIFICATE OF REGISTRATION

This is to certify that

## Promega Biosciences, LLC

277 Granada Drive, San Luis Obispo, California 93401, USA

operates a

## Quality Management System

which complies with the requirements of

## ISO 9001:2015

for the following scope of certification

**Design and Manufacture of Specialty Chemicals.**

Certificate No.: CERT-0135851  
File No.: 1045467  
Issue Date: July 7, 2020

Original Certification Date: April 5, 2005  
Certification Effective Date: August 8, 2020  
Certificate Expiry Date: August 7, 2023

Heather Mahon  
Global Head of Technical Services  
SAI Global Assurance



ISO 9001



Registered by:  
QMI-SAI Canada Limited (SAI Global), 20 Carlson Court, Suite 200, Toronto, Ontario M9W 7K6 Canada. This registration is subject to the SAI Global Terms and Conditions for Certification. While all due care and skill was exercised in carrying out this assessment, SAI Global accepts responsibility only for proven negligence. This certificate remains the property of SAI Global and must be returned to them upon request.  
To verify that this certificate is current, please refer to the SAI Global On-Line Certification Register:  
[https://www.saiglobal.com/en-us/assurance/auditing\\_and\\_certification/certification\\_registry/](https://www.saiglobal.com/en-us/assurance/auditing_and_certification/certification_registry/)

 **SAI GLOBAL**  
INFORM. INSPIRE. IMPROVE.