



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-0104462
File No.: 1045467
Issue Date: August 9, 2017

Original Certification Date: April 5, 2005
Certification Effective Date: August 9, 2017
Certificate Expiry Date: August 8, 2020

Nicole Grantham
General Manager SAI Global Certification Services



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: www.qmi-saiglobal.com/qmi_companies/

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