



Certificate

The SQS herewith attests that the organisation named below has a management system that meets the requirements of the normative base mentioned.



Promega AG
Wallisellenstrasse 55
8600 Dübendorf
Switzerland

Scope

**Supply of biological/chemical reagents
and scientific instruments**

Normative base

ISO 9001:2015

Quality Management System

Reg. no. 14538

Validity 17.11.2019–16.11.2022
Issue 17.11.2019


A. Grisard, President SQS


F. Müller, CEO SQS



sqs.ch



Swiss Association for Quality
and Management Systems (SQS)
Bernstrasse 103, 3052 Zollikofen, Switzerland

