



# Certificate

SQS herewith certifies that the company named below has a management system which meets the requirements of the standard specified below.



**Promega AG**  
**Wallisellenstrasse 55**  
**8600 Dübendorf**  
**Switzerland**

Scope of certification

**Whole Company**

Field of activity

**Supplier of biological reagents, scientific instruments  
and disposables**

Normative base

**ISO 9001:2015**

**Quality Management System**

Scope(s) 12

Validity 17.11.2016 – 16.11.2019  
Issue 17.11.2016

Reg. no. 14538

X. Edelmann, President SQS

R. Glauser, CEO SQS



sqs.ch



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

