



# 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  
Page 1 of 1

Validity 05.12.2025 – 16.11.2028  
Issue 05.12.2025

4020\_2 / Januar2024\_Version 3.1



*A. Grisard*  
A. Grisard, President SQS

*F. Müller*  
F. Müller, CEO

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

