



THE INTERNATIONAL CERTIFICATION NETWORK

CERTIFICATE

SQS has issued an IQNet recognized certificate that the organization:

Sartorius Stedim Switzerland AG

Ringstrasse 24a

8317 Tagelswangen

Switzerland

has implemented and maintains a

Management System

for the following scope(s):

18

which fulfills the requirements of the following standard(s):

ISO 9001:2015 Quality Management System

Issued on: 2019-04-12

Expires on: 2022-04-11

This attestation is directly linked to the IQNet Partner's original certificate and shall not be used as a stand-alone document

Registration Number: CH-45289



Alex Stoichitoiu
President of IQNet

Felix Müller
CEO SQS



IQNet Partners*:

AENOR Spain AFNOR Certification France APCER Portugal CCC Cyprus CISQ Italy
CQC China CQM China CQS Czech Republic Cro Cert Croatia DQS Holding GmbH Germany FCAV Brazil
FONDONORMA Venezuela ICONTEC Colombia Inspecta Sertifointi Oy Finland INTECO Costa Rica
IRAM Argentina JQA Japan KFQ Korea MIRTEC Greece MSZT Hungary Nemko AS Norway NSAI Ireland
NYCE-SIGE México PCBC Poland Quality Austria Austria RR Russia SII Israel SIQ Slovenia
SIRIM QAS International Malaysia SQS Switzerland SRAC Romania TEST St Petersburg Russia TSE Turkey YUQS Serbia
IQNet is represented in the USA by: AFNOR Certification, CISQ, DQS Holding GmbH and NSAI Inc.

* The list of IQNet partners is valid at the time of issue of this certificate. Updated information is available under www.iqnet-certification.com



Certificate

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



sartorius

Sartorius Stedim Switzerland AG
Ringstrasse 24a
8317 Tagelswangen
Switzerland

Scope of certification

Whole Company

Field of activity

**Development and production
of products for biotechnology**

Normative base

ISO 9001:2015

Quality Management System

Scope(s) 18

Validity 12.04.2019 – 11.04.2022
Issue 12.04.2019

Reg. no. 45289

X. Edelmann, President SQS

F. Müller, CEO SQS



sqs.ch



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





Zertifikat

Die SQS bescheinigt hiermit, dass nachstehend genanntes Unternehmen über ein Managementsystem verfügt, das den Anforderungen der aufgeführten normativen Grundlage entspricht.



sartorius

Sartorius Stedim Switzerland AG

Ringstrasse 24a

8317 Tagelswangen

Schweiz

Geltungsbereich

Ganzes Unternehmen

Tätigkeitsgebiet

**Entwicklung und Produktion
von Produkten für die Biotechnologie**

Normative Grundlage

ISO 9001:2015

Qualitätsmanagementsystem

Scope(s) 18

Gültigkeit 12.04.2019 – 11.04.2022
Ausgabe 12.04.2019

Reg.-Nr. 45289

X. Edelmann, Präsident SQS

F. Müller, CEO SQS



sqs.ch



Schweizerische Vereinigung für
Qualitäts- und Management-Systeme SQS
Bernstrasse 103, 3052 Zollikofen, Schweiz

