

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Tecan Schweiz AG
Seestraße 103
8708 Männedorf
Switzerland

Holds Certificate No:

FM 666054

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.

For and on behalf of BSI:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2014-05-28

Latest Revision Date: 2022-06-13

Effective Date: 2022-06-19

Expiry Date: 2025-06-18

Page: 1 of 2



...making excellence a habit.™

Certificate No: **FM 666054**

Location	Registered Activities
Tecan Schweiz AG Seestraße 103 8708 Männedorf Switzerland	The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.
Tecan Schweiz AG Emil Staub Str. 1 8708 Männedorf Switzerland	The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.
Tecan Schweiz AG Gewerbstrasse 8 8606 Nänikon Switzerland	The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.



Original Registration Date: 2014-05-28

Latest Revision Date: 2022-06-13

Effective Date: 2022-06-19

Expiry Date: 2025-06-18

Page: 2 of 2

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.
An electronic certificate can be authenticated [online](#).
Printed copies can be validated at www.bsigroup.com/ClientDirectory

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 345 080 9000

BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.