

# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Tecan Trading AG  
Seestraße 103  
Männedorf  
8708  
Switzerland

Holds Certificate Number:

FM 666071

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

**Sales, distribution, installation and service of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.**

For and on behalf of BSI:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2014-05-28

Latest Revision Date: 2020-05-15

Effective Date: 2020-05-28

Expiry Date: 2023-05-27

Page: 1 of 2



...making excellence a habit.™

Certificate No: FM 666071

Location

Tecan Trading AG  
Seestraße 103  
Männedorf  
8708  
Switzerland

Tecan Trading AG  
Emil Staub Str. 1  
Männedorf  
8708  
Switzerland

Registered Activities

Sales, distribution, installation and service of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.

Sales, distribution, installation and service of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.



Original Registration Date: 2014-05-28

Latest Revision Date: 2020-05-15

Effective Date: 2020-05-28

Expiry Date: 2023-05-27

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](http://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.  
A Member of the BSI Group of Companies.