

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016 & EN ISO 13485:2016

This is to certify that:

Tecan Schweiz AG

Seestraße 103
8708 Männedorf
Switzerland

Holds Certificate Number:

MD 666058

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 & EN ISO 13485:2016 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:



Stewart Brain, Head of Compliance & Risk - Medical Devices

Original Registration Date: 2017-02-20

Latest Revision Date: 2018-06-06

Effective Date: 2017-05-01

Expiry Date: 2019-02-28

Page: 1 of 2



003

...making excellence a habit.™

Certificate No: MD 666058

Location

Tecan Schweiz AG
Seestraße 103
8708 Männedorf
Switzerland

Tecan Schweiz AG
Emil Staub Str. 1
8708 Männedorf
Switzerland

Registered Activities

Design and Development of robotic sample processors and automated solutions for in-vitro diagnostics.

Manufacturing of robotic sample processors and automated solutions for in-vitro diagnostics.



Original Registration Date: 2017-02-20

Latest Revision Date: 2018-06-06

Effective Date: 2017-05-01

Expiry Date: 2019-02-28

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.
A Member of the BSI Group of Companies.