

Certificate of Registration

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

This is to certify that:

Tecan Trading AG
Seestraße 103
Männedorf
8708
Switzerland

Holds Certificate Number:

MD 666074

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

Sales, distribution, installation and service 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: 2017-02-16

Latest Revision Date: 2018-06-06

Effective Date: 2017-05-01

Expiry Date: 2020-04-30

Page: 1 of 2



003

...making excellence a habit.™

Certificate No: **MD 666074**

Location

Registered Activities

Tecan Trading AG
Seestraße 103
Männedorf
8708
Switzerland

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

Tecan Trading AG
Emil Staub Str. 1
Männedorf
8708
Switzerland

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



Original Registration Date: 2017-02-16

Effective Date: 2017-05-01

Latest Revision Date: 2018-06-06

Expiry Date: 2020-04-30

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.