

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 Operations and Compliance Director

Original Registration Date: 28/05/2014

Latest Revision Date: 03/04/2017

Effective Date: 28/05/2017

Expiry Date: 27/05/2020



003

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: 28/05/2014

Latest Revision Date: 03/04/2017

Effective Date: 28/05/2017

Expiry Date: 27/05/2020