

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Tecan Austria GmbH
Untersbergstraße 1a
5082 Grödig
Austria

Holds Certificate Number:

FM 675530

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

Design and Development, Production, Sales, Distribution, Installation and Service of Devices for General Laboratory Use and Devices for In-vitro Diagnostic (including Accessories)

For and on behalf of BSI:



Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2015-01-12

Latest Revision Date: 2018-01-18

Effective Date: 2017-12-20

Expiry Date: 2020-12-19



003

Page: 1 of 1

...making excellence a habit.™