

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Tecan Austria GmbH
Untersbergstraße 1a
5082 Grödig
Austria

Holds Certificate Number:

MD 675529

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

The design, development, manufacture, distribution, service of instruments and accessories for in-vitro diagnostics.



For and on behalf of BSI:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2017-07-12

Latest Revision Date: 2020-07-07

Effective Date: 2020-12-20

Expiry Date: 2023-06-03

Page: 1 of 1



003

...making excellence a habit.™