

# 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:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2017-07-12

Latest Revision Date: 2023-05-23

Effective Date: 2023-06-04

Expiry Date: 2026-06-03

Page: 1 of 1



...making excellence a habit.™