

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: 2020-06-18

Effective Date: 2020-06-18

Expiry Date: 2023-06-03

Page: 1 of 1



003

...making excellence a habit.™