

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016 & EN ISO 13485:2016

This is to certify that:

Tecan Software Competence Center GmbH
Peter Sander Strasse 41a
Mainz-Kastel
55252
Germany

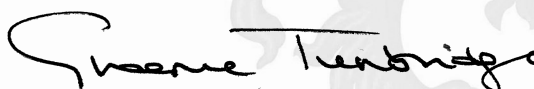
Holds Certificate Number:

MD 723201

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

Design, development and maintenance of software for robotic sample processors and automated solutions for in-vitro diagnostics.

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2020-09-14

Latest Revision Date: 2023-08-24

Effective Date: 2023-09-14

Expiry Date: 2026-09-13

Page: 1 of 1



...making excellence a habit.™