

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Tecan US Inc.
9401 Globe Center Drive
Suite 140
Morrisville
North Carolina
27560
USA

Holds Certificate Number:

MD 759104

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

Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in- vitro diagnostic applications. Sales and distribution of reagents and consumables.

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2022-09-19

Latest Revision Date: 2022-09-19

Effective Date: 2022-09-19

Expiry Date: 2025-09-18

Page: 1 of 1



...making excellence a habit.™