

Certificate of Registration

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

This is to certify that:

Tecan SP, Inc.
14180 Live Oak Ave
Baldwin Park
CA 91706
USA

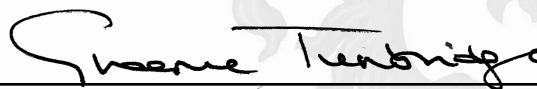
Holds Certificate Number:

MD 694547

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

The design and manufacture of instruments and consumables, and the installation and services of instruments for solid phase extraction for medical and in- vitro diagnostic use.

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Global Regulatory & Quality

Original Registration Date: 2018-12-06

Latest Revision Date: 2024-11-07

Effective Date: 2024-12-06

Expiry Date: 2027-12-05

Page: 1 of 1



...making excellence a habit.™