

# 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 for solid phase extraction for medical and in-vitro diagnostic use.

For and on behalf of BSI:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2018-12-06

Latest Revision Date: 2021-12-01

Effective Date: 2021-12-06

Expiry Date: 2024-12-05

Page: 1 of 1



...making excellence a habit.™