

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 used in life science.

For and on behalf of BSI:



Stewart Brain, Head of Compliance & Risk - Medical Devices

Original Registration Date: 2018-12-06

Latest Revision Date: 2018-12-06

Effective Date: 2018-12-06

Expiry Date: 2021-12-05



003

Page: 1 of 1

...making excellence a habit.™