

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

Tecan Systems Inc.

2450 Zanker Road
San Jose
California
95131-1126
USA

Holds Certificate Number:

MD 670834

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

The design, development, manufacture and service of pump components and robotic liquid handling systems for life science applications.

For and on behalf of BSI:



Stewart Brain, Head of Compliance & Risk - Medical Devices

Original Registration Date: 2017-06-02

Latest Revision Date: 2017-08-04

Effective Date: 2017-08-31

Expiry Date: 2020-08-30

Page: 1 of 1



003

...making excellence a habit.™