

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Tecan Systems Inc.
2450 Zanker Road
San Jose
California
95131-1126
USA

Holds Certificate No:

FM 670835

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

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

This certificate is traceable to this company's original certificate number QS5 14 09 54551 006 issued by TUV SUD.

For and on behalf of BSI:



Carlos Pitanga, SVP, System Certification and Compliance

Original Registration Date: 2014-09-03

Latest Revision Date: 2017-07-25

Effective Date: 2017-09-03

Expiry Date: 2020-09-02

Page: 1 of 1



...making excellence a habit.™