

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.

For and on behalf of BSI:

Carlos Pitanga, Chief Operating Officer Assurance – Americas

Original Registration Date: 2014-09-03

Latest Revision Date: 2020-11-25

Effective Date: 2020-09-03

Expiry Date: 2023-09-02

Page: 1 of 1



...making excellence a habit.™