

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

Tecan Systems Inc.
18635 Sutter Boulevard
Morgan Hill
California
95037
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 provision of design and development services as well as servicing of pump components, robotic liquid handling systems and integrated instruments for life science applications.

For and on behalf of BSI:

Carlos Pitanga, Managing Director Assurance, Americas

Original Registration Date: 2014-09-03

Latest Revision Date: 2023-08-17

Effective Date: 2023-09-03

Expiry Date: 2026-09-02

Page: 1 of 2



...making excellence a habit.™

Certificate No: **FM 670835**

Location

Registered Activities

Tecan Systems Inc.
18635 Sutter Boulevard
Morgan Hill
California
95037
USA

The provision of design and development services as well as servicing of pump components, robotic liquid handling systems and integrated instruments for IVD (in-vitro diagnostic) and medical device applications.

Tecan Systems Inc.
18735 Madrone Parkway
Morgan Hill
California
California
95037
USA

Servicing of pump components, robotic liquid handling systems and integrated instruments for IVD (in-vitro diagnostic) and medical device applications.



Original Registration Date: 2014-09-03

Effective Date: 2023-09-03

Latest Revision Date: 2023-08-17

Expiry Date: 2026-09-02