

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:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2017-06-02

Latest Revision Date: 2022-09-16

Effective Date: 2020-08-31

Expiry Date: 2023-08-30



Page: 1 of 2

...making excellence a habit.™

Certificate No: **MD 670834**

Location

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

Registered Activities

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

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

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



Original Registration Date: 2017-06-02

Latest Revision Date: 2022-09-16

Effective Date: 2020-08-31

Expiry Date: 2023-08-30

Page: 2 of 2

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.
An electronic certificate can be authenticated [online](#).
Printed copies can be validated at www.bsigroup.com/ClientDirectory

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 345 080 9000
BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.
A Member of the BSI Group of Companies.