

# Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that:

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

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2017-06-02

Latest Revision Date: 2023-08-15

Effective Date: 2023-08-31

Expiry Date: 2026-08-30



Page: 1 of 2

...making excellence a habit.™

Certificate No: **MD 670834**

Location

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

Registered Activities

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: 2017-06-02

Latest Revision Date: 2023-08-15

Effective Date: 2023-08-31

Expiry Date: 2026-08-30

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](http://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.