

# 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 Number:

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, Managing Director Assurance, Americas

Original Registration Date: 2014-09-03

Latest Revision Date: 2022-11-20

Effective Date: 2020-09-03

Expiry Date: 2023-09-02

Page: 1 of 2



...making excellence a habit.™

Certificate No: FM 670835

Location

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

Registered Activities

The design, development, 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

Design and development activities



Original Registration Date: 2014-09-03

Latest Revision Date: 2022-11-20

Effective Date: 2020-09-03

Expiry Date: 2023-09-02

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.