

# Certificate of Registration

## QUALITY MANAGEMENT SYSTEM - ISO 13485:2016

This is to certify that: Tecan Schweiz AG  
Seestraße 103  
8708 Männedorf  
Switzerland

Facility ID Number: F002717

Holds Certificate No: **MDSAP 705120**

Statement of Conformity: The company listed on this certificate has been audited to and found to conform with the following criteria: ISO 13485:2016 and Australia - Therapeutic Goods (Medical Devices) Regulations, 2002, Schedule 3 Part 1 (excluding Part 1.6) - Full Quality Assurance Procedure; Brasil - RDC ANVISA n. 16/2013, RDC ANVISA n. 23/2012, RDC ANVISA n. 67/2009; Canada - Medical Devices Regulations - Part 1 - SOR 98/282; Japan - MHLW Ministerial Ordinance 169, Article 4 to Article 68, PMD Act; USA - 21 CFR 820, 21 CFR 803, 21 CFR 806, 21 CFR 807 - Subparts A to D

Design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.

For and on behalf of BSI:

Gary E Slack, Senior Vice President - Medical Devices

Original Registration Date: 2019-06-19

Effective Date: 2021-06-10

Expiry Date: 2022-06-18



BSI Group America Inc. is an MDSAP authorized auditing organization

Page: 1 of 2

...making excellence a habit.™

Certificate No: **MDSAP 705120**

Location	Registered Activities
Tecan Schweiz AG Seestraße 103 8708 Männedorf Switzerland Facility ID Number: F002717	Design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.
Tecan Schweiz AG Emil Staub Str. 1 8708 Männedorf Switzerland Facility ID Number: F002717	Design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.
Tecan Schweiz AG Gewerbstrasse 8 8606 Nänikon Switzerland Facility ID Number: F002717	Incoming inspection and warehousing of components and sub assemblies for robotic sample processors and automated solutions for in-vitro diagnostics.



Original Registration Date: 2019-06-19

Effective Date: 2021-06-10

Expiry Date: 2022-06-18

Page: 2 of 2

This certificate remains the property of BSI and shall be returned immediately upon request.  
An electronic certificate can be authenticated [online](http://www.bsigroup.com/ClientDirectory). Printed copies can be validated at [www.bsigroup.com/ClientDirectory](http://www.bsigroup.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.

Americas Headquarters: BSI Group America Inc., 12950 Worldgate Drive, Suite 800, Herndon, VA 20170-6007 USA  
A Member of the BSI Group of Companies.