

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016 & EN ISO 13485:2016

This is to certify that:

Tecan Schweiz AG

Seestraße 103
8708 Männedorf
Switzerland

Holds Certificate Number:

MD 806640

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 & EN ISO 13485:2016 for the following scope:

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

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Global Regulatory & Quality

Original Registration Date: 2024-05-21

Latest Revision Date: 2024-05-21

Effective Date: 2024-05-21

Expiry Date: 2025-06-18

Page: 1 of 2



...making excellence a habit.™

Certificate No: MD 806640

Location	Registered Activities
Tecan Schweiz AG Seestraße 103 8708 Männedorf Switzerland	The 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	The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.
Tecan Schweiz AG Gewerbstrasse 8 8606 Nänikon Switzerland	The design, development and manufacture of robotic sample processors and automated solutions for in-vitro diagnostics.



Original Registration Date: 2024-05-21

Latest Revision Date: 2024-05-21

Effective Date: 2024-05-21

Expiry Date: 2025-06-18

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 Group The Netherlands B.V., John M. Keynesplein 9, 1066 EP Amsterdam, The Netherlands | Tel: +31 20 3460 780

BSI Group The Netherlands B.V. is registered in The Netherlands under number 33264284 | A Member of the BSI Group Holdings B.V.