

Certificate of Registration

ENVIRONMENTAL MANAGEMENT SYSTEM - ISO 14001:2015

This is to certify that:

Tecan Trading AG
Seestrasse 103
Männedorf
8708
Switzerland

Holds Certificate No:

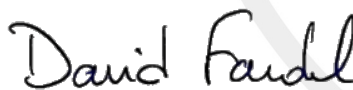
EMS 785658

and operates an Environmental Management System which complies with the requirements of ISO 14001:2015 for the following scope:

The design, development, manufacture, warehouse, sales, distribution, installation and robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.

Activities are related to the IAF Sector 19.

For and on behalf of BSI:



David Fardel, Country Manager, Assurance - Continental Europe

Original Registration Date: 2023-09-29

Latest Revision Date: 2023-09-29

Effective Date: 2023-09-29

Expiry Date: 2026-09-28



Page: 1 of 2

...making excellence a habit.™

Certificate No: **EMS 785658**

Location	Registered Activities
Tecan Trading AG Seestrasse 103 Männedorf 8708 Switzerland	The design, development, sales, distribution, installation and service of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.
Tecan CH Seestrasse 103 Männedorf 8708 Switzerland	The manufacture of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.
Tecan CH Gewerbstrasse 8 Nänikon 8606 Switzerland	The warehouse of robotic sample processors and automated solutions for in-vitro diagnostics and life science service solutions.

Original Registration Date: 2023-09-29

Latest Revision Date: 2023-09-29

Effective Date: 2023-09-29

Expiry Date: 2026-09-28

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 contacts: BSI Group Italia srl, via Fara 35, Milano, Italy, phone: +39 02 6679091
A Member of the BSI Group of Companies.