

Certificate of Registration

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

This is to certify that:

Tecan Deutschland GmbH
Werner-von-Siemens-Straße 23
Crailsheim
74564
Germany

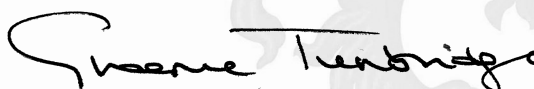
Holds Certificate Number:

MD 676093

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

Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2017-09-25

Latest Revision Date: 2024-01-18

Effective Date: 2024-01-31

Expiry Date: 2027-01-30

Page: 1 of 3



...making excellence a habit.™

Certificate No: **MD 676093**

Location	Registered Activities
Tecan Deutschland GmbH Werner-von-Siemens-Straße 23 Crailsheim 74564 Germany	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan France S.A.S.U Tour Swiss Life 1 Boulevard Marius Vivier Merle Lyon 69003 France	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan Benelux bvba Schaliënhoevedreef 20A Mechelen 2800 Belgium	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan Sales Austria GmbH Untersbergstrasse 1a Grödig 5082 Austria	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Sales Office Tecan Nordic AB Sveavägen 159 1tr Stockholm 11346 Sweden	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan Ibérica Instrumentación S.L. C/Lepanto 151 Bajos Barcelona 08013 Spain	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan Italia S. r. l. Via Brescia 39 Cernusco sul Naviglio MI 20063 Italy	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.

Original Registration Date: 2017-09-25

Latest Revision Date: 2024-01-18

Effective Date: 2024-01-31

Expiry Date: 2027-01-30

Page: 2 of 3

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

Certificate No: **MD 676093**

Location	Registered Activities
Tecan Sales Switzerland Seestrasse 103 Männedorf 8708 Switzerland	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan UK Ltd. Theale Court 11-13 High Street Theale RG7 5AH United Kingdom	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan Australia Pty Ltd 21/3 Westside Avenue Port Melbourne Victoria 3207 Australia	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.
Tecan (Shanghai) Laboratory Equipment Co.,Ltd. 12F, 1529 Guozhan Road Pudong New Area Shanghai 200122 China	Sales, distribution, installation and service of robotic sample processors, detection devices and automated laboratory solutions for in-vitro diagnostic applications.

Original Registration Date: 2017-09-25

Latest Revision Date: 2024-01-18

Effective Date: 2024-01-31

Expiry Date: 2027-01-30

Page: 3 of 3

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, 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.