



CERTIFICATE



This is to certify that

Tecan Trading AG

Seestrasse, 103
8708 Männedorf
Switzerland

with the organizational units/sites as listed in the annex

has implemented and maintains an **Information Security Management System**.

Scope:

The information Security System of Tecan Trading AG campus to ensure security and privacy for the provision of development, production and distribution of automated workflow solutions for laboratories in the life sciences sector.

With reference to Statement of Applicability (SOA): Ver 2.0, 2026-06-03.

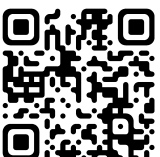
Through an audit, documented in a report, it was verified that the management system fulfills the requirements of the following standard:

ISO / IEC 27001 : 2022

Certificate registration no.	31633375 ISMS22
Date of original certification	2026-08-26
Date of certification	2026-08-26
Valid until	2029-08-25



DQS IS A MEMBER OF



DQS Inc.

David Tellez
Managing Director

Accredited Body: DQS Inc., 1500 McConnor Parkway, Suite 400, Schaumburg, IL 60173 USA
The validity of the certification can only be verified by the QR-code.



Annex to certificate
Registration no.: 31633375 ISMS22

Tecan Trading AG

Seestrasse, 103
8708 Männedorf
Switzerland

Location

10002924
Tecan CDMO Solutions MH, Corp.
18735 Madrone Parkway
Morgan Hill, CA 95037
United States of America

31628792
Tecan CDMO Solutions MH, Corp.
18715 Madrone Parkway
Morgan Hill, CA 95037
United States of America

31637454
Tecan Switzerland AG
Emil-Staub Strasse 1
8708 Männedorf
Switzerland

31637455
Tecan Switzerland AG
Gewerbestrasse 8
8606 Nänikon
Switzerland

Scope

The information Security System of Tecan CDMO campus to ensure security and privacy for the production and distribution of automated workflow solutions for laboratories in the life sciences sector. With reference to the SOA Ver 2.0, 2026-06-03.

The information Security System of Tecan CDMO campus to ensure security and privacy for the production and distribution of automated workflow solutions for laboratories in the life sciences sector. With reference to the SOA Ver 2.0, 2026-06-03.

The information Security System of Tecan Trading AG campus to ensure security and privacy for the provision of development, production and distribution of automated workflow solutions for laboratories in the life sciences sector. With reference to the SOA Ver 2.0, 2026-06-03.

The information Security System of Tecan Trading AG campus to ensure security and privacy for the provision of development, production and distribution of automated workflow solutions for laboratories in the life sciences sector. With reference to the SOA Ver 2.0, 2026-06-03.