



CERTIFICATE

This is to certify that

Tecan CDMO Solutions MH, Corp.

18735 Madrone Parkway
Morgan Hill, CA 95037
United States of America

with the organizational units/sites as listed in the annex

has implemented and maintains a **Quality Management System**.

Scope:

Design, manufacture and service of non-implantable electronic components and box-build assemblies.

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

ISO 9001 : 2015

Certificate registration no.	10002924 QM15
Date of original certification	1995-04-25
Date of revision	2025-09-08
Date of certification	2024-05-22
Valid until	2027-05-26



DQS Inc.

David Tellez
Managing Director

DQS IS A MEMBER OF



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.: 10002924 QM15

Tecan CDMO Solutions MH, Corp.

18735 Madrone Parkway
Morgan Hill, CA 95037
United States of America

Location

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

Scope

Product Realization, customer related processes,
Production and service provision.