



CERTIFICATE



This is to certify that

Tecan CDMO Solutions MH, Corporation

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

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

Scope:

Design, manufacture and service of non-implantable electronic components, assemblies and finished medical devices.

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

ISO 13485 : 2016

Certificate registration no.	10002924 MP2016
Date of original certification	2006-04-22
Date of revision	2024-12-23
Date of certification	2024-05-22
Valid until	2027-05-26



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.