



# CERTIFICATE



This is to certify that

## Paramit Corporation (A Tecan Group Company)

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	2024-05-23
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.