



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 an **Information Security Management System**.

Scope:

The Information Security Management System (ISMS) of Paramit Corporation, a fully owned company of Tecan Group, is to ensure information Security and Privacy for the provision of services covering the confidentiality, integrity and availability of the operations and customers information maintained in our information processing systems.

With reference to Statement of Applicability (SOA): 30703TMt02 V1.3, 2024-05-20

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.	10002924 ISMS22
Date of original certification	2024-07-09
Date of certification	2024-07-09
Valid until	2027-07-08



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.