



America

CERTIFICATE

No. QS2 066559 0003 Rev. 01

Certificate Holder:

Tecan Genomics
900 Chesapeake Drive, Ground Floor
Redwood City CA 94063
USA

Certification Mark:



Scope of Certificate:

Design, Development, Manufacture and Distribution of Nucleic Acid Amplification Technology for use in Scientific Research and Medical Diagnostic Fields

Standard:

ISO 13485:2016

The Certification Body of TÜV SÜD America Inc. certifies that the company mentioned above has established and is maintaining a quality management system that meets the requirements of the listed standards.

Report No.:

72156549

Effective Date:

2020-08-30

Expiry Date:

2023-04-18

Page 1 of 1

Date of Issue: 2020-09-26

(Tina Israel)
Manager, US Certification Body,
Medical and Health Services