



Certificate

No. Q8 070179 0007 Rev. 00

Holder of Certificate: Tecan Japan Co., Ltd.
Kawasaki Tech Center
580-16, Horikawacho, Saiwai-ku
Kawasaki-shi
Kanagawa
212-0013 JAPAN



Facility(ies): Tecan Japan Co., Ltd.
Kawasaki Tech Center, 580-16, Horikawacho,
Saiwai-ku, Kawasaki-shi, Kanagawa, 212-0013 JAPAN

Certification Mark:



Scope of Certificate: Sales, distribution, installation, and servicing of
in-vitro diagnostic instruments for clinical
diagnosis

Applied Standard(s): EN ISO 13485:2016
Medical devices - Quality management systems -
Requirements for regulatory purposes
(ISO 13485:2016)
DIN EN ISO 13485:2016

The Certification Body of TÜV SÜD Product Service GmbH certifies that the company mentioned above has established and is maintaining a quality management system, which meets the requirements of the listed standard(s). See also notes overleaf.

Report No.: JAQ235034211

Valid from: 2018-10-01
Valid until: 2021-09-30

Date, 2018-08-28

Stefan Preiß

ZERTIFIKAT ◆ CERTIFICATE ◆ 認證證書 ◆ CERTIFICADO ◆ CERTIFICAT