

Declaration of Software Verification

FLUENTCONTROL 3.6

FluentControl 3.6 software has been designed, verified and validated according to the Tecan Product Development Process. Software Verification is performed to ensure that the Design Output meets the Design Input Requirements. Design Validation is performed to ensure that the product meets the User Needs and Intended Use. Typical process documents of the Product Development Process include but are not limited to Verification Reports, Design Validation Reports, Risk Management, as well as Manuals and Help files. Master Media for software distribution is scanned to be malware-free.

Tecan performs IQ and OQ on instrument installation. PQ is the customer's responsibility and is specific to their Fluent's application.

Our commitment to continuous improvement is certified by our registration to ISO 9001:2015 and ISO 13485:2016 quality system standards. Tecan's software development complies with IEC 62304 – medical device software – software life cycle processes. To view our certificates online, please scan the QR code below.

Rachel Begheyn

Product Manager Software, T-CH

DocuSigned by:



Signer Name: Rachel Begheyn
Signing Reason: I am the author of this document
Signing Time: 2024-07-05 | 4:15:42 PM CEST

7E4EDC397849491AA71464AA4CFDE30E

Asmaa Edres-Lohaus

Software Project Manager, T-SCC

DocuSigned by:



Signer Name: Asmaa Edres-Lohaus
Signing Reason: I have reviewed this document
Signing Time: 2024-07-05 | 4:23:15 PM CEST

A9ED6834BB2A4728B05D5BD3DEE21266

Johann Israel

Director QA DACH, T-CH

DocuSigned by:



Signer Name: Johann Israel
Signing Reason: I approve this document
Signing Time: 2024-07-05 | 4:19:33 PM CEST

1B7C02D2F58F42FD89226D4DA71CF4B3

