

EU DECLARATION OF CONFORMITY

This declaration of conformity is issued under the sole responsibility of the manufacturer:

Tecan Schweiz AG, Seestrasse 103, CH-8708 Männedorf

for the product:

Freedom EVO 75

Part No.	Model
30025019	FREEDOM EVO 75
30025869	FREEDOM EVO 75
Software:	Freedom EVOware® Standard (for 30025019) Freedom EVOware® Plus (for 30025869)
Options:	Liquid Handling Arm™ Robotic Manipulator Arm™

is in conformity with the provisions of the following European Directive(s) when installed in accordance with the installation instructions contained in the product documentation:

Directive 2006/42/EC

on machinery

Directive 2014/30/EU

relating to electromagnetic compatibility

Directive 2011/65/EU

on the restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS 2) including Commission Delegated Directive (EU) 2015/863 (RoHS3) amending Annex II to Directive 2011/65/EU

and that the standards referenced below were taken in consideration:

EN 61010-1:2010+A1:2019

Safety requirements for electrical equipment for measurement, control and laboratory use - Part 1: General requirements

EN 61010-2-010:2020

Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-010: Particular requirements for laboratory equipment for the heating of material

EN 61010-2-081:2020

Safety requirements for electrical equipment for measurement, control and laboratory use - Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes

EN 61326-1:2021

Electrical equipment for measurement, control and laboratory use - EMC requirements - Part 1: General Requirements

EN 62304: 2006+A1:2015

Medical Device software – Software life cycle processes

EN 62366-1: 2015

Medical Device software – Application of usability engineering to medical devices

EN ISO 15223-1: 2021Medical devices. Symbols to be used with medical device labels, labelling and information to be supplied.
General requirements**EN ISO 14971: 2012**

Medical devices – Application of risk management to medical devices

EN ISO 12100: 2010

Safety of machinery - General principles for design - Risk assessment and risk reduction

EN IEC 63000: 2018

Technical documentation for the assessment of electrical and electronic products with respect to the restriction of hazardous substances

Tecan Schweiz AG maintains a quality system certified to the following standards:

EN ISO 9001: 2015

Quality management systems – Requirements

EN ISO 13485: 2016

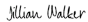
Medical devices – quality Management Systems – Requirements for regulatory purposes

Tecan Schweiz AG, Seestrasse 103, CH-8708 Männedorf 2025-02-17

Signed by:

Signer Name: Johann Israel
Signing Reason: I approve this document
Signing Time: 2025-02-17 | 6:22:57 PM CET
1B7C02D2F58F42FD89226D4DA71CF4B3

Johann Israel
Director QA DACH

Signed by:

Signer Name: Jillian Walker
Signing Reason: I approve this document
Signing Time: 2025-02-18 | 8:30:40 AM CET
9CBF24ABC93B424D81F6BBF9FE61FE0B

Jillian Walker
Associate Director RA Instrumentation

Certificate Of Completion

Envelope Id: 5C52005C-C97C-4073-A651-8742334F1CA8	Status: Completed
Subject: Complete with Docusign: 396926.pdf EU DoC EVO 75	
Source Envelope:	
Document Pages: 2	Signatures: 2
Certificate Pages: 2	Initials: 0
AutoNav: Enabled	Envelope Originator:
Envelopeld Stamping: Disabled	Adriana Franco
Time Zone: (UTC+01:00) Amsterdam, Berlin, Bern, Rome, Stockholm, Vienna	Seestrasse 103
	Männedorf, ZH 8708
	Adriana.Franco@Tecan.com
	IP Address: 91.212.29.30

Record Tracking

Status: Original	Holder: Adriana Franco	Location: DocuSign
2/17/2025 5:54:02 PM	Adriana.Franco@Tecan.com	

Signer Events

Signer Events	Signature	Timestamp
Jillian Walker		Sent: 2/17/2025 5:58:13 PM
Jillian.walker@tecan.com		Viewed: 2/18/2025 8:30:20 AM
Associate Director Regulatory Affairs		Signed: 2/18/2025 8:30:45 AM
Instrumentation		
Tecan Schweiz AG	Signature Adoption: Pre-selected Style	
Security Level: Email, Account Authentication (Required)	Signature ID: 9CBF24AB-C93B-424D-81F6-BBF9FE61FE0B	
	Using IP Address: 91.212.29.29	
	With Signing Authentication via Docusign password	
	With Signing Reasons (on each tab): I approve this document	

Electronic Record and Signature Disclosure:

Not Offered via Docusign

Johann Israel		Sent: 2/17/2025 5:58:13 PM
Johann.Israel@tecan.com		Viewed: 2/17/2025 6:22:35 PM
Director QA Tecan.		Signed: 2/17/2025 6:23:04 PM
Tecan Schweiz AG	Signature Adoption: Pre-selected Style	
Security Level: Email, Account Authentication (Required)	Signature ID: 1B7C02D2-F58F-42FD-8922-6D4DA71CF4B3	
	Using IP Address: 89.26.17.169	
	Signed using mobile	
	With Signing Authentication via Docusign password	
	With Signing Reasons (on each tab): I approve this document	

Electronic Record and Signature Disclosure:

Not Offered via Docusign

In Person Signer Events	Signature	Timestamp
Editor Delivery Events	Status	Timestamp
Agent Delivery Events	Status	Timestamp
Intermediary Delivery Events	Status	Timestamp
Certified Delivery Events	Status	Timestamp
Carbon Copy Events	Status	Timestamp

Witness Events	Signature	Timestamp
Notary Events	Signature	Timestamp
Envelope Summary Events	Status	Timestamps
Envelope Sent	Hashed/Encrypted	2/17/2025 5:58:13 PM
Certified Delivered	Security Checked	2/17/2025 6:22:35 PM
Signing Complete	Security Checked	2/17/2025 6:23:04 PM
Completed	Security Checked	2/18/2025 8:30:45 AM
Payment Events	Status	Timestamps