

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:

Fluent[®]

Part No.	Model
30042011	INSTRUMENT FLUENT 480 BASE UNIT
30042021	INSTRUMENT FLUENT 780 BASE UNIT
30042031	INSTRUMENT FLUENT 1080 BASE UNIT
Software:	FluentControl™
Options:	Robotic Gripper Arm™ (RGA) Robotic Gripper Arm™ Long Z (RGA-Z) Multiple Channel Arm™ 384/96 (MCA384) Multiple Channel Arm™ (MCA) 96 with optional Gripper Flexible Channel Arm™ (FCA) Air Flexible Channel Arm™ (Air FCA) Air Flexible Channel Arm™ (Air FCA) Multisense Fluent ID™ Fluent Stacker™ with/ without Barcode reader Fluent Carousel™ Monitored Incubator Option™ (MIO-2) Te-Shake™ Te-VacS™ Tube Rotator™ Frida Reader™ ResolveX® i300

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 2014/30/EU

relating to electromagnetic compatibility

Directive 2006/42/EC

on machinery

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 IEC 61010-2-051: 2021+A11:2021

Safety requirements for electrical equipment for measurement, control and laboratory use - Part 2-051: Particular requirements for laboratory equipment for mixing and stirring

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 60825-1: 2014+A11:2021

Safety of laser products – Part 1: Equipment classification and requirements

EN 62304: 2006+A1:2015

Medical Device software – Software life cycle processes

EN 62366-1: 2015+A1:2020

Medical Device software – Application of usability engineering to medical devices

EN ISO 15223-1: 2021

Medical devices. Symbols to be used with medical device labels, labelling and information to be supplied. General requirements

EN ISO 18113-1: 2011

In vitro diagnostic medical devices - Information supplied by the manufacturer (labelling) - Part 1: Terms, definitions and general requirements

EN ISO 18113-3: 2011

In vitro diagnostic medical devices -- Information supplied by the manufacturer (labelling) -- Part 3: In vitro diagnostic instruments for professional use

EN ISO 14971: 2019+A11:2021

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

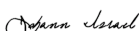
Quality management systems – Requirements

EN ISO 13485: 2016

Medical devices – quality Management Systems – Requirements for regulatory purposes

Tecan Schweiz AG, Seestrasse 103, CH-8708 Männedorf 2024-08-12

Signed by:



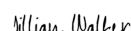

Signer Name: Johann Israel

Signing Reason: I approve this document

Signing Time: 2024-08-12 | 1:25:56 PM CEST

1B7C02D2F58F42FD89226D4DA71CF4B3

Signed by:




Signer Name: Jillian Walker



Signing Reason: I approve this document

Signing Time: 2024-08-12 | 1:31:06 PM CEST

9CBF24ABC93B424D81F6BBF9FE61FE0B

Johann Israel
Director QA

Jillian Walker
Associate Director RA Instrumentation

Certificate Of Completion		
Envelope Id: 5FAEEDF1FD26472DA28E3A0863079E5D	Status: Completed	
Subject: Complete with DocuSign: 398421 V12.2_draft.docx		
Source Envelope:		
Document Pages: 2	Signatures: 2	Envelope Originator:
Certificate Pages: 2	Initials: 0	Kurt Erwin Hurab
AutoNav: Enabled		Seestrasse 103
Envelopeld Stamping: Disabled		Männedorf, ZH 8708
Time Zone: (UTC+01:00) Amsterdam, Berlin, Bern, Rome, Stockholm, Vienna		KurtErwin.Hurab@tecan.com
		IP Address: 81.223.7.138
Record Tracking		
Status: Original	Holder: Kurt Erwin Hurab	Location: DocuSign
2024-08-12 13:21	KurtErwin.Hurab@tecan.com	
Signer Events	Signature	Timestamp
Jillian Walker		Sent: 2024-08-12 13:23
Jillian.walker@tecan.com		Viewed: 2024-08-12 13:30
Associate Director Regulatory Affairs		Signed: 2024-08-12 13:31
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: 62.2.160.130	
	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		
Johann Israel		Sent: 2024-08-12 13:23
Johann.Israel@tecan.com		Viewed: 2024-08-12 13:25
Director QA Tecan.		Signed: 2024-08-12 13:26
Tecan Schweiz AG	Signature Adoption: Pre-selected Style	
Security Level: Email, Account Authentication (Required)	Signature ID: 1B7C02D2-F58F-42FD-8922-6D4DA71CF4B3	
	Using IP Address: 62.2.160.130	
	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	2024-08-12 13:23
Certified Delivered	Security Checked	2024-08-12 13:25
Signing Complete	Security Checked	2024-08-12 13:26
Completed	Security Checked	2024-08-12 13:31

Payment Events	Status	Timestamps
----------------	--------	------------