

UK 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 UK Regulations when installed in accordance with the installation instructions contained in the product documentation:

Electromagnetic Compatibility Regulations 2016
The Supply of Machinery (Safety) Regulations 2008
The Restriction of the Use of Certain Hazardous Substances in
Electrical and Electronic Equipment Regulations 2012

and that the standards referenced below were taken in consideration:

BS EN 61010-1: 2010+A1:2019

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

BS 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

BS 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

BS 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

BS EN 61326-1: 2021

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

BS EN 60825-1: 2014+A11:2021

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

BS EN 62304: 2006+A1:2015

Medical Device software – Software life cycle processes

BS EN 62366-1: 2015+A1:2020

Medical Device software – Application of usability engineering to medical devices

BS EN ISO 15223-1: 2021

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

BS EN ISO 18113-1: 2011

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

BS 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

BS EN ISO 14971: 2019 + A11:2021

Medical devices – Application of risk management to medical devices

BS EN ISO 12100: 2010

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

BS EN IEC 63000: 2018

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

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

BS EN ISO 9001: 2015

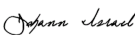
Quality management systems – Requirements

BS EN ISO 13485: 2016

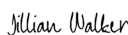
Medical devices – quality Management Systems – Requirements for regulatory purposes

Person authorized to compile the technical file: Tecan UK Ltd, Theale Court, 11 – 13 High Street, Theale, Reading, RG7 5AH, United Kingdom

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 | 1:26:39 PM CEST
1B7C02D2F58F42FD89226D4DA71CF4B3

Johann Israel
Director QA

Signed by:

 Signer Name: Jillian Walker
Signing Reason: I approve this document
Signing Time: 2024-08-12 | 1:32:44 PM CEST
9CBF24ABC93B424D81F6BBF9FE61FE0B

Jillian Walker
Associate Director RA Instrumentation

Certificate Of Completion		
Envelope Id: 36CE9C3C74AA42EB9A4D34C9BFAE2CB9		Status: Completed
Subject: Complete with DocuSign: 402020 V3.2_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 13:24	KurtErwin.Hurab@tecan.com	
Signer Events	Signature	Timestamp
Jillian Walker		Sent: 2024-08-12 13:25
Jillian.walker@tecan.com		Viewed: 2024-08-12 13:31
Associate Director Regulatory Affairs		Signed: 2024-08-12 13:32
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 13:25
Johann.Israel@tecan.com		Viewed: 2024-08-12 13:26
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 13:25
Certified Delivered	Security Checked	2024-08-12 13:26
Signing Complete	Security Checked	2024-08-12 13:26
Completed	Security Checked	2024-08-12 13:32

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