

EU DECLARATION OF CONFORMITY

This declaration of conformity is issued under the sole responsibility of the manufacturer:
Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria

for the product:

SPARK

Part No.(or Cat. No.)	Model	Configuration
30086376	SPARK	Injector, Heater and Magnetic Stirrer

Valid from serial number: **2405000001**

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 60825-1: 2014+A11:2021

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

EN 61326-1: 2021

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

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 Austria GmbH maintains a quality system certified to the following standards:

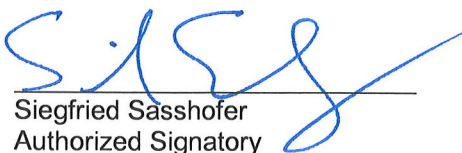
EN ISO 9001: 2015

Quality management systems – Requirements

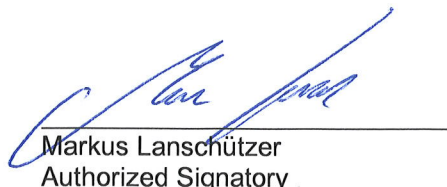
EN ISO 13485: 2016

Medical devices – Quality management systems – Requirements for regulatory purposes

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria, 2024-05-14



Siegfried Sasshofer
Authorized Signatory



Markus Lanschützer
Authorized Signatory

UK DECLARATION OF CONFORMITY

This declaration of conformity is issued under the sole responsibility of the manufacturer:
Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria

for the product:

SPARK

Part No.(or Cat. No.)	Model	Configuration
30086376	SPARK	Injector, Heater and Magnetic Stirrer

Valid from serial number: **2405000001**

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

The 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 60825-1: 2014 + A11:2021

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

BS EN 61326-1: 2021

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

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 Austria GmbH 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, Court, 11 – 13 High Street, Theale, Reading, RG7 5AH, United Kingdom

Tecan Austria GmbH Untersbergstr. 1A A-5082 Grödig, Austria, 2024-05-14



Siegfried Sasshofer
Authorized Signatory



Markus Lanschützer
Authorized Signatory