



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Flow Injection System, Model FIAS 100

This is to certify that this PerkinElmer product was tested and verified to be in conformance with all applicable quality requirements, including specifications, drawings, calibration, preservation, packing, marking requirements and part identification.

Declaration of EMC, Safety, and RoHS Compliance

This PerkinElmer product conforms to the regulations stipulated in the CE Mark requirements for the EMC Directive (2014/30/EU), the Low Voltage Directive (2014/35/EU), and the RoHS 2 Directive (2011/65/EU):

EN 55011:2009 + A1:2010, Group 1, Class A, EMC -- RF Characteristics of ISM Equipment
EN 61326-1:2013, EMC -- Requirements for Electrical Equipment for Laboratory Use
EN 61000-4-2:2009, EMC -- Electrostatic Discharge Requirements
EN 61000-4-3:2006 + A1:2008 + A2:2010, EMC -- Radiated Electromagnetic Field Requirements
EN 61000-4-4:2004 + A1:2010, EMC -- Electrical Fast Transient/Burst Requirements
EN 61000-4-5:2006, EMC -- Surge Immunity Requirements
EN 61000-4-6:2009, EMC -- Conducted Disturbances (induced by RF fields) Requirements
EN 61000-4-8:2010, EMC -- Power Frequency Magnetic Field Requirements
EN 61000-4-11:2004, EMC -- Voltage Dips, Short Interruptions, Voltage Variations Requirements
EN 61000-3-2:2006 + A1:2009 + A2:2009, EMC -- Harmonic Current Emissions
EN 61000-3-3:2008, EMC -- Voltage Fluctuations and Flicker
EN 61326-2-2:2013, EMC -- Particular requirements for portable and low voltage equipment
EN 61010-1:2010, Safety Requirements for Electrical Equipment for Laboratory Use
EN 50581:2012, Technical documentation for the assessment of electrical and electronic products with respect to the RoHS



CAN/CSA C22.2 No. 61010-1-12, Safety Requirements for Electrical Equipment for Laboratory Use
UL 61010-1, 3rd edition, Safety Requirements for Electrical Equipment for Laboratory Use
CAN/CSA C22.2 No. 61010-2-081:15, Particular Requirements for Automatic & Semi-automatic Laboratory Equipment for Analysis & Other Purposes
UL 61010-2-081, 2nd edition, Particular Requirements for Automatic and Semi-automatic Laboratory Equipment for Analysis and Other Purposes
ICES-003 Issue 6, Class A, Radiated and Conducted Emissions
FCC Part 15, Class A, Radiated and Conducted Emissions

AS/NZS CISPR 11:2011
Korean Radio Waves Act, Article 58-2, Clause 3

NOTE: The operation of certain types of equipment (e.g., signal generators) may be subject to given restrictions. Please refer to the appropriate information in the respective user documentation.

Declaration of System Validation

The product was found to meet its functional and performance specification prior to shipment. To support this declaration, the following Engineering, Assembly and Test documents are held by PerkinElmer and are available for reference upon request in justified cases and to an appropriate extent:

The Product Description	The System Design Documentation
The Functional Specification	The Source Code Documentation
The User Interface Definition	The Evaluation Documentation

NOTE: PerkinElmer will maintain possession of all documents and controls their reproduction, including parts of them.

The existence of these documents and the procedures used in their production are formal requirements of the PerkinElmer Quality Management System. The integrity of the PerkinElmer Quality Management System is routinely audited and has been certified to ISO 9001 since 1992.

This declaration of conformity is issued under the sole responsibility of PerkinElmer.

Signed for and on behalf of:

Alan Mears
Compliance Engineer

22 April 2020

09936583G.doc

PerkinElmer, 710 Bridgeport Avenue, Shelton, CT 06484 USA
[in the EU: PerkinElmer, Chalfont Road, Buckinghamshire, Seer Green, HP9 2FX, United Kingdom]
An ISO 9001 Company