



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Flexar Solvent Manager

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: 2012, 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 61010-1:2012, Safety Requirements for Electrical Equipment for Laboratory Use
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
ICES-003, Class A, Radiated and Conducted Emissions
AS/NZS CISPR 11:2011
Korean Radio Waves Act, Article 58-2, Clause 3
FCC Part 15, Class A, Radiated and Conducted Emissions



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 Functional Specification
The User Interface Definition

The System Design Documentation
The Source Code Documentation
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:

Rodney Wright
Compliance Engineer

09 September, 2017

09936844B.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