



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

### Flexar LC Pumps

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 61326-1:2012, EMC -- Requirements for Electrical Equipment for Laboratory Use  
EN 61000-4-2:2008, 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 61000-4-5:2005, EMC -- Surge Immunity Requirements  
EN 61000-4-6:2008, EMC -- Conducted Disturbances (induced by RF fields) Requirements  
EN 61000-4-8:2009, EMC -- Power Frequency Magnetic Field Requirements  
EN 61000-4-11:2009, 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, 3<sup>rd</sup> 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 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

17 July 2019