



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

### Atomic Absorption Spectrometer, Models PinAAcle 900F, 900H, 900T, and 900Z

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 61326-2-1:2013, EMC -- Particular requirements for sensitive test and measurement equipment for EMC unprotected applications  
EN 61000-3-2:2006 + A1:2009 + A2:2009, EMC -- Harmonic Current Emissions  
EN 61000-3-3:2008, EMC -- Voltage Fluctuations and Flicker  
EN 61010-1:2010, Safety Requirements for Electrical Equipment for Laboratory Use  
EN 61010-2-010:2014, Safety Requirements for Laboratory Equipment for the Heating of Materials  
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, 3<sup>rd</sup> edition, Safety Requirements for Electrical Equipment for Laboratory Use  
CAN/CSA C22.2 No. 61010-2-010:15, Safety Requirements for Laboratory Equipment for the Heating of Materials  
UL 61010-2-010, 3<sup>rd</sup> edition, Safety Requirements for Laboratory Equipment for the Heating of Materials  
CAN/CSA C22.2 No. 61010-2-061-15, Safety Requirements for Laboratory Spectrometers with Thermal Atomization and Ionization  
UL 61010-2-061, 3<sup>rd</sup> edition, Safety Requirements for Laboratory Spectrometers with Thermal Atomization and Ionization  
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, 2<sup>nd</sup> 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  
Law on Metrology of the People's Republic of China (Pattern Approval; Model PinAAcle 900T only)



*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 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:

Alan Mears  
Compliance Engineer

21 April 2020