



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Gas Mixing Device, Model GMD 8000

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 61000-4-5:2015, EMC -- Surge Immunity Requirements
 - EN 61000-4-6:2014, 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 61000-3-2:2014, EMC -- Harmonic Current Emissions
- EN 61000-3-3:2013, EMC -- Voltage Fluctuations and Flicker
- 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
- ICES-003, Class B, Radiated and Conducted Emissions
- FCC Part 15, Class B, 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

11 June 2019