



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Gas Chromatograph, Model Clarus 590

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), the Machinery Directive (2006/42/EC), 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 61000-4-11:2004, 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 61010-2-010:2014, Safety Requirements for Laboratory Equipment for the Heating of Materials
 - EN 61010-2-081:2015, Particular Requirements for Automatic and Semi-automatic Laboratory Equipment for Analysis and Other Purposes
 - 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 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:


Rodney Wright
Compliance Engineer

05 March 2018



An ISO 9001 Company