

Declaration of Conformity

PerkinElmer, Inc.
68 Elm Street
Hopkinton, MA 01748
USA

Declares that the following product:

Type of Equipment:

Electrical Laboratory Equipment

Model Name and Part Number(s):

Model Name

Quantum GX2 Instrument

Part Numbers

CLS149276

Has been manufactured using the following standards and specifications to comply with:

Directives
2014/35/EU (Low Voltage)
2014/30/EU (Electro Magnetic Compatability)
2011/65/EU (RoHS)

Electrical Safety
IEC/EN 61010-1: 2010 Third Edition
CAN/CSA C22.2 No. 61010-1:2012
UL 61010-1:2012

Restriction of the use of certain Hazardous Substances (RoHS)
EN 50581: 2012

Electromagnetic compatability
EN 61326-1: 2013
EN 55011:2009 +A1:2010, Group 1, Class A
KN 11
KN 61000-6-1

National standards and technical specification:

None

We have an internal production control system that ensures compliance between the manufactured products and the Technical Construction File. PerkinElmer, Inc. holds the TCF. Refer to the Bill of Materials for the above part number(s) for Technical Construction File information.

Declaration of Conformity

As manufacturer we declare, under our sole responsibility that the equipment follows the provisions of the standards and specifications stated above.

Signature: Lisa Allen
Date: 13-APR-2018
Title of signatory: Sr. Quality Manager