

Declaration of Conformity

PerkinElmer
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

Part Numbers

IVIS Spectrum

124262, 124374, 124375

IVIS Spectrum BL

CLS137325, CLS137326, CLS137327

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

Directives and Amendments:

2014/35/EU	Low Voltage Directive
2014/30/EU	EMC Directive
768/2008/EC	CE Marking Directive
2011/65/EU	RoHS directive

Harmonized standards used:

EN 61010-1: 2010	Electrical Safety
EN 55011:2009A/A1:2010, Group 1 Class A	EMI for Industrial, Scientific and Medical Equipment
EN 61326-1: 2013	EMC Requirements / Electrical Equipment for Measurement, Control and Laboratory Use
EN 55011:2009/A1:2010, Group 1, Class A	EMI for Industrial, Scientific and Medical Equipment

National standards and technical specification:

Australia & New Zealand	ACMA AS/NZS CISPR 11:2009/A1:2010, Group 1 Class A	EMC Requirements / Industrial scientific and medical (ISM) radio-frequency equipment
Canada	CAN/CSA C22.2 No. 61010-1:2012	Electrical Safety
	ICES-003, Issue 6, Class A	EMC for Information Technology Equipment (Including Digital Apparatus)
USA	CFR Title 47 FCC Part 15 Subpart B, Class A	EMC for Radio Frequency Devices – Unintentional Radiators
	UL 61010-1:2012	Electrical Safety for Electrical Equipment For Measurement, Control, and Laboratory Use

Declaration of Conformity

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

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

Signature:

-disa all

Date:

03-NOV-2017

Title of signatory:

Sr. Quality Manager