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
EN 33011.2009/A1.2010, Group 1, Class A	Medical Equipment

National standards and technical specification:

	1 with statute as and technical specification.		
Australia &	ACMA AS/NZS CISPR	EMC Requirements / Industrial	
New	11:2009/A1:2010, Group 1	scientific and medical (ISM) radio-	
Zealand	Class A	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	EMC for Radio Frequency Devices –	
	Subpart B, Class A	Unintentional Radiators	
		Electrical Safety for Electrical	
	UL 61010-1:2012	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: Title of signatory:

atory: Sr. Quality Manager