



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

IVIS Lumina X5 Imaging Systems, Part Number CLS148590

Type of Equipment: Electrical Laboratory Equipment

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 61000-4-4:2004 + A1:2010, 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 61000-4-8:2010, EMC -- Power Frequency Magnetic Field Requirements
EN 61000-4-11:2004, EMC -- Voltage Dips, Short Interruptions, Voltage Variations Requirements
EN 61000-3-2:2014, Class A, EMC -- Harmonic Current Emissions
EN 61000-3-3:2013, Voltage Fluctuations and Flicker
EN 61010-1:2010, Safety Requirements for Electrical Equipment for Laboratory Use

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, Issue 6, Class A, Radiated and Conducted Emissions
FCC Part 15, Subpart B, Class A, Radiated and Conducted Emissions

ACMA AS/NZS CISPR 11:2011, Group 1, Class A

Korean Radio Waves Act, Article 58-2, Clause 3



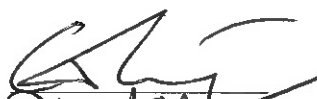
This declaration of conformity is issued under the sole responsibility of PerkinElmer.

Signed for and on behalf of:

Signature:

Title of signatory:

Date:


Dir. of QA
05/15/2018

PerkinElmer, 68 Elm Street, Hopkinton, MA, USA

P/N CLS150271 Rev A