

# *RoHS Certificate of Compliance*



PerkinElmer.  
68 Elm Street  
Hopkinton, Massachusetts 01748  
(508) 435-9500  
FAX (508) 435-3439

## ***CLS137325 IVIS Instrument, Spectrum BL, 120V***

### **Restriction of the use of certain Hazardous Substances (RoHS)**

All items in this assembly do not exceed the maximum concentration values (MCVs) in EU Directive 2011/65/EC as stated in Annex II of that directive, and therefore, this assembly is RoHS Compliant.

This means that each of the homogenous materials within this product does not exceed the following:

- Lead (0.1 %)
- Mercury (0.1 %)
- Cadmium (0.01 %)
- Hexavalent chromium (0.1 %)
- Polybrominated biphenyls (PBB) (0.1 %)
- Polybrominated diphenyl ethers (PBDE) (0.1 %)