

RoHS Certificate of Compliance



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

CLS137326 IVIS Instrument, Spectrum BL, 100V

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 %)