

RoHS Certificate of Compliance



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

124374 ***IVIS Instrument, Spectrum, 100V, Andor Camera***

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 %)