RoHS Certificate of Compliance



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

124375 IVIS Instrument, Spectrum, 220V, 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 %)

CLS145008 Rev. A Page 1 of 1