## RoHS Certificate of Compliance



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

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

The RoHS Directive 2011/65/EU restricts the use of certain hazardous substances in electrical and electronic equipment.

Items that support the production of PerkinElmer equipment but are not included as part of the actual instrument are considered RoHS compliant. The following are included in this category:

- Procedures
- Documents
- Tools
- Software
- Graphic Images
- Reference Items
- Quality Control Materials

CLS145007 Rev. A Page 1 of 1