

Declaration of REACH Compliance

Product Description: Infrared Spectroscopy
 Model Number: Spotlight 150 and Spotlight 200
 Assembly Number:

PerkinElmer, Inc. declares that the products listed above do contain Substances of Very High Concern (SVHCs) as defined in the Candidate List of the REACH Regulation EC 1907/2006. REACH SVHCs which are not found on the Authorization list are not restricted but must be declared in accordance with REACH Article 33. Based on the information available the products contain the following SVHC(s)

Substance	CAS	EC
1,2-dimethoxyethane; ethylene glycol dimethyl ether (EGDME)	110-71-4	203-794-9
1,3,5-Tris(oxiran-2-ylmethyl)-1,3,5-triazinane-2,4,6-trione (TGIC)	2451-62-9	219-514-3
2-(2H-benzotriazol-2-yl)-4,6-ditertpentylphenol (UV-328)	25973-55-1	247-384-8
4,4'-isopropylidenediphenol	80-05-7	201-245-8
Benzene-1,2,4-tricarboxylic acid 1,2 anhydride	552-30-7	209-008-0
Bis(hexachlorocyclopentadieno)cyclooctane	13560-89-9	236-948-9
Boric acid	10043-35-3	233-139-2
Boric acid	11113-50-1	234-343-4
Cobalt dichloride	7646-79-9	231-589-4
Cyclohexane-1,2-dicarboxylic anhydride	85-42-7	201-604-9
Diboron trioxide	1303-86-2	215-125-8
Lead	7439-92-1	231-100-4
Lead monoxide (lead oxide)	1317-36-8	215-267-0
Octamethylcyclotetrasiloxane	556-67-2	209-136-7

This declaration is based on PerkinElmer, Inc. understanding of the requirements of the REACH Regulation and knowledge of the materials that go into its products. PerkinElmer, Inc. bases its knowledge on information provided by third-party suppliers and makes no representation or warranty as to the accuracy of such information. PerkinElmer, Inc. continues to take steps to obtain accurate information from suppliers but has not conducted destructive testing or chemical analysis on incoming materials to verify material composition.

Date Issued:

10/25/19
 PerkinElmer, Inc.
 710 Bridgeport Avenue
 Shelton, CT 06484-4794 United States of America

Authorized Signature:

