

Declaration of REACH Compliance

Product Description: Liquid Chromatography
Model Number: QSight 110 and QSight 120
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-Methyl-2-pyrrolidone	872-50-4	212-828-1
Benzene-1,2,4-tricarboxylic acid 1,2 anhydride	552-30-7	209-008-0
Bis (2-ethylhexyl) phthalate (DEHP)	117-81-7	204-211-0
Diboron trioxide	1303-86-2	215-125-8
Hexahydro-1-methylphthalic anhydride	48122-14-1	256-356-4
Hexahydro-3-methylphthalic anhydride	57110-29-9	260-566-1
Hexahydro-4-methylphthalic anhydride	19438-60-9	243-072-0
Hexahydromethylphthalic anhydride	25550-51-0	247-094-1
Lead	7439-92-1	231-100-4
Lead monoxide (lead oxide)	1317-36-8	215-267-0
Lead titanium zirconium oxide	12626-81-2	235-727-4
Refractory ceramic fibers	142844-00-6	

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:

