

Declaration of REACH Compliance

Product Description: UV/VIS Spectroscopy

Model Number: Lambda 465 UV-VIS

Assembly Number:

PerkinElmer, Inc. declares that, to its knowledge as of the date of this document, the products listed above conform to the requirements of the EU REACH Regulation EC 1907/2006, and to the best of our knowledge we have determined that none of the Substances of Very High Concern (SVHCs) listed in the July 16, 2019 Candidate List are present over the 0.1% by weight threshold in any article contained in the product.

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:

