



2020 Disclosure
United Kingdom Modern Slavery Act 2015
Bio-Rad Laboratories, Inc.

This disclosure by Bio-Rad Laboratories, Inc. is made on behalf of its subsidiaries that provide goods or services in the United Kingdom.

Bio-Rad Laboratories, Inc. is a United States-based corporation that manufactures and supplies the life science research, healthcare, analytical chemistry and other markets with a broad range of products and systems used to separate complex chemical and biological materials and to identify, analyze and purify their components. Bio-Rad Laboratories, Inc. operates in two industry segments designated as Life Science and Clinical Diagnostics, and both segments operate worldwide. In 2020, Bio-Rad Laboratories, Inc. had three subsidiaries operating in the United Kingdom that supplied goods and services -- Bio-Rad Laboratories Limited, Bio-Rad AbD Serotec Ltd, and Bio-Rad Services UK Limited. The Global Supply Chain organization of Bio-Rad Laboratories, Inc. operates worldwide and has responsibility for the manufacturing, procurement and distribution functions that support the three subsidiaries in the United Kingdom.

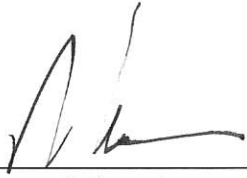
The policy of Bio-Rad Laboratories, Inc. in relation to slavery and human trafficking, implemented through its Global Supply Chain organization, is that all labor employed in the production of its products must be voluntary as stated in its Supplier Code of Conduct. Bio-Rad Laboratories, Inc. requires that its suppliers do not engage in or support trafficking in human beings. Slave, child, underage, forced, bonded, or indentured labor will not be tolerated by Bio-Rad Laboratories, Inc. Upon the request of Bio-Rad Laboratories, Inc., its suppliers shall certify that they have implemented procedures to ensure compliance with laws on slavery and human trafficking.

The Global Supply Chain organization of Bio-Rad Laboratories, Inc. evaluates the risk of slavery and human trafficking in the supply chain and also has procedures to mitigate potential risks. Use of a third-party compliance tool has been initiated to further assist the Global Supply Chain organization in identifying potential supplier risks related to modern-day slavery, employee rights, and employee health and safety practices. The Global Supply Chain organization also conducts periodic audits of its principal suppliers.

Bio-Rad Laboratories, Inc. makes available to its workforce members, officers, directors, business partners and other third parties (except where prohibited by law) a "helpline" through which concerns or violations of policy or unethical conduct – including those associated with slavery, human trafficking or other labor abuses – may be reported and further investigated by Bio-Rad Laboratories, Inc. Such disclosure through the helpline is confidential and anonymous. The availability of this communication channel, along with appropriate instructions, is provided

in Bio-Rad Laboratories, Inc.'s Supplier Code of Conduct and its Code of Business Ethics and Conduct.

Bio-Rad Laboratories, Inc.'s Code of Business Ethics and Conduct states that purchasers engage only those suppliers who comply with all laws regarding human trafficking and slavery. All of Bio-Rad Laboratories, Inc.'s employees and employees of its subsidiaries undergo periodic training on the Code of Business Ethics and Conduct. The Global Supply Chain organization of Bio-Rad Laboratories, Inc. provides periodic training to employees responsible for procurement of materials and services and has undertaken to familiarize certain personnel with identifying slavery and human trafficking issues and mitigating such risks.



Date: May 7, 2021

Norman Schwartz

Chairman of the Board, President and Chief Executive Officer, Bio-Rad Laboratories, Inc.
Director, Bio-Rad Laboratories Limited
Director, Bio-Rad Services UK Limited
Director, Bio-Rad AbD Serotec Ltd.