

SARTORIUS

Responsibility

UK Modern
Slavery Statement

Simplifying Progress

This statement provides information about supply chain transparency at Sartorius and has been published in accordance with the U.K. Modern Slavery Act 2015 (the “Act”). It sets out the steps that Sartorius and its affiliates, including The Automation Partnership (Cambridge) Limited to which the Act applies, have taken during the financial year ending 31 December 2020 and plan to continue taking, to prevent modern slavery in its businesses and direct supply chains.

Unless otherwise specified, all references to Sartorius herein are references to the above-named Sartorius entities and to Sartorius globally to the extent it is subject to the Act.

Modern slavery encompasses slavery, servitude, human trafficking and forced labour. Sartorius has a zero tolerance approach to any form of modern slavery. Sartorius is committed to acting ethically and with integrity and transparency in all business dealings and to putting effective systems and controls in place to safeguard against any form of modern slavery taking place within the business or its supply chain.

About Sartorius:

Sartorius is a leading international pharmaceutical and laboratory equipment supplier. With innovative products and services, Sartorius is helping its customers across the entire globe to implement complex and quality-critical bio-manufacturing and laboratory processes reliably and economically.

The Group companies are united under the roof of Sartorius AG, which is listed on the Frankfurt Stock Exchange and holds the majority stake in Sartorius Stedim Biotech S.A. quoted on the Paris Stock Exchange, this subgroup is comprised mainly of the Bioprocess Solutions Division to which The Automation Partnership (Cambridge) Limited belongs.

Sartorius Code of Conduct for Suppliers

Sartorius not only supplies high-quality products, but also does so in a manner that reflects the corporation's mission and commitment to conducting its business activities in compliance with the prevailing laws in a humane, socially responsible, and sustainable way.

Sartorius expects that all of its business partners, and particularly its suppliers, behave likewise. The Sartorius Code of Conduct and the Sartorius Anti-Corruption Code set forth a defined set of values and principles to which Sartorius categorically subscribes. The present Supplier Code of Conduct has been written to promote continual implementation of these codicils by setting forth a set of incontrovertible minimum standards. Sartorius obliges its suppliers and their employees, representatives, agents and subcontractors (hereinafter referred to as Suppliers) to observe and uphold these standards when conducting business activities.

Compliance with International Social and Environmental Standards

Sartorius suppliers are required to uphold international agreements and guidelines, including the Universal Declaration of Human Rights, the conventions of the International Labour Organization (ILO) and the United Nations Global Compact.

The relevant standards are available on the United Nation's website: www.un.org.

Suppliers are required, among others, to ensure:

- Environmental Protection
- Health and Safety in the Workplace
- Equal Opportunity
- Prohibition of Forced Labour
- Prohibition of Child Labour
- Remuneration to meet legal minimum wage and/or industry standards

Should Sartorius find any evidence that these codes have been violated by a Supplier, it will undertake to prosecute any breaches by suitable means.

Compliance and Training

With its compliance management systems that is valid worldwide, Sartorius ensures that the members of its individual boards, executives and employees comply with all legal regulations and codes, and perform their activities in accordance with the company's internal rules and guidelines. Targeted training and awareness raising prevent any misconduct, as well as economic damage and loss of image.

Sartorius makes every effort to ensure optimal risk management by using a combination of approaches: a preventive compliance approach designed to proactively stop any potential breaches before they occur and a repressive compliance approach intended to continuously monitor compliance with the company's rules. These processes are closely intermeshed, creating a standardized compliance management system that aims to offer the best possible protection against potential violations of rules and regulations. Sartorius has developed a Code of Conduct as a preventive component of its compliance management system and has committed to an Anti- Corruption Code. An internal system is available for reporting any suspicious circumstances involving potential compliance violations.

This statement has been reviewed and endorsed by the Board of Directors of The Automation Partnership (Cambridge) Limited.



Scott Fegan

Director

8th June 2021