



Sartorius Code of Conduct for suppliers

March 2013

Preamble

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 Anticorruption Code of Conduct 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. We oblige our employees, representatives, agents and subcontractors (hereinafter referred to as Suppliers) to observe and uphold these standards when conducting business activities.

It is the responsibility of Suppliers to instruct their employees, representatives, agents and subcontractors accordingly. Should Sartorius find any evidence that these codes have been violated, it will undertake to prosecute any breaches by suitable means.

1 Compliance with International Social and Environmental Standards

Supplier shall uphold international agreements and guidelines, including the Universal Declaration of Human Rights, the conventions of the International Labor Organization (ILO) and the United Nations Global Compact.

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

Environmental Protection

Supplier shall help protect the environment by developing, manufacturing and distributing its products in a resource-saving manner. The ecological effects of our business activities shall be regularly monitored and assessed.

Health and Safety in the Workplace

Supplier shall be committed to offering all its employees a safe and healthy working environment and to continually improving this environment.

Equal Opportunity

In hiring and employing human resources, Supplier shall not discriminate against, disadvantage, harass or exclude anyone based on gender, ethnic origin, life philosophy, race, religion, age, disability, appearance, sexual preferences and identity, origin or political position.

Prohibition of Forced Labor

Supplier is prohibited from using any forced, bonded or otherwise indentured labor.

Prohibition of Child Labor

Child labor and any form of exploitation of children by Supplier are prohibited. Supplier shall respect the special need to protect young employees.

Remuneration

The remuneration paid for regular working hours, overtime and compensation for hours worked in excess of contract or regular work schedules shall meet the legal minimum wage and/or industry standards. Remuneration may not be withheld illegally or as a means of punishment. Remuneration shall be paid in accordance with the laws in force.

The Right to Collective Bargaining and Freedom of Association

Supplier shall respect the right of all employees to form and join unions and worker representative groups. Supplier's employees who are members of a union or of a worker representative group shall neither be treated preferentially nor be disadvantaged.

2 General Rules of Conduct

Compliance with Laws and Statutes

We hold the compliance with laws and statutes to be self-evident. Accordingly, Sartorius expects its Suppliers always to adhere to the terms and conditions of the legal framework of the countries in which they operate.

Fair Competition

In line with fair competition, suppliers shall observe and uphold all applicable anti-trust laws, which include but are not limited to a ban on price fixing, market share and capacity agreements, the division of regional markets or customers and controlled pricing.

3 Dealing with Conflicts of Interest

Suppliers shall not influence Sartorius employees in their awarding of contracts by offering perks.