

SARTORIUS

Responsibility

Sartorius'
Environmental Position

Simplifying Progress

Contents

1	Environmental protection in the corporate strategy	3
2	Fundamental principles of our environmental position	3
3	Core topics in our environmental position	4

Document information

Applicability	
Scope	Sartorius Group
Confidentiality	Public
Ownership	
Policy Owner	Katharina Tillmanns
Board Area	Group Strategy, Human Resources, Corporate Research, Legal & Compliance, Communications
Board Sponsor	Joachim Kreuzburg
Approved by	Joachim Kreuzburg
Document information	
Release	February 2019
Review	Bi-annually

1 Environmental protection in the corporate strategy

Sartorius has offices and production facilities in more than 110 countries on six continents. It is an integral part of our corporate strategy that we pursue our business activities in a resource-efficient and thus environmentally responsible manner.

We are a signatory of the United Nations Global Compact and support its Sustainable Development Goals (SDGs). Our internal guidelines and processes are oriented toward the objective of limiting our impact on the environment in an economically and socially appropriate way.

Sartorius' Environmental Position is binding on all the companies and sites of the Sartorius Group. It has been communicated to all our employees via the intranet, and supports us in anchoring efficiency and environmental awareness in our daily business. The management of Sartorius is accountable for the implementation of the environmental position. The Executive Board sets goals on the basis of the fundamental principles and core topics in the company's position paper. Both global and local measures are carried out to realize all the environmentally specific objectives. For this reason, certified environmental management systems will be put in place successively at all of the Group's manufacturing facilities. As a matter of principle, even sites that have not been certified are operated according to the same high standards.

2 Fundamental principles of our environmental position

Compliance

Regular audits ensure our compliance with statutory regulations and requirements. If a country's laws prescribe no requirements – or standards that are less stringent than our internal guidelines – action plans are developed to close the gap step by step.

Transparency, dialog and ongoing improvement

We want to create transparency as regards our environmental impact and promote a common understanding of environmental protection in our company. This is why we pursue an open dialog with our employees and business partners as well as with state institutions. The aim of this dialog is the ongoing improvement of our operational environmental protection in harmony with the expectations of our stakeholders.

Employees

We keep our employees informed about environmental protection topics and encourage them to act responsibly. In addition, we train them on environmental topics relevant to their individual tasks.

Inclusion of the value creation chain

We work in tandem with customers and suppliers to arrive at a common understanding of environmental protection. We expect our suppliers to implement these principles, which are also set down in our Code of Conduct for Suppliers.

High environmental protection standards

It is our aspiration to keep the environmental impact of our activities as low as possible. The use of natural resources and the generation of waste should rise disproportionately lower than sales. We drive forward technological progress to improve the environmental footprint of our products, and we use the opportunities we have to help solve regional and global environmental problems.

Products

Our products must meet the safety and quality requirements of our customers in the biopharmaceutical industry. At the same time, we aim for a high degree of environmental compatibility in our products.

3 Core topics in our environmental position

Sartorius has defined its core topics with respect to environmental protection. These are topics of particular importance to our daily business and our stakeholders. We have separate internal guidelines and instructions for these core topics.

Single-use products

Single-use products enable manufacturing processes that are both safe and flexible – and thus economical. In addition, production with single-use products frequently consumes fewer resources than traditional production using stainless steel equipment. Single-use products can thus significantly reduce the environmental impact of biopharmaceutical processes. Sartorius rigorously considers the environmental footprint of its single-use products in all their life cycle phases as early as the product development stage: The materials used are carefully selected to keep their impact on humans and the environment as low as possible. The plastics consist of recyclable standard polymers and contain few additives. Whenever technically possible, the materials meet Pharmacopeia requirements. All materials can be safely burnt and used as energy.

Water and hazardous materials

Our manufacturing processes for membranes and membrane products are optimized for minimal consumption of rinsing water. Organic solvents are processed and largely recycled. We comply with German and international regulations in our use and handling of hazardous materials, and we always pursue the goal of forgoing hazardous materials when we develop new products.

Locations and building technology

We invest continuously in new plants and plant expansions to accommodate our constant growth. In constructing new facilities worldwide, we follow recognized standards for sustainable building as guidance.

Transportation and packaging

Our company-wide commitment to long-term environmental protection also includes our logistics. As early as the product development phase, we choose the sizes and materials of our packaging as per environmental protection criteria. In particular, we use materials made of renewable raw materials (such as cardboard) as well as films and plastics that are 100% recyclable. We also apply ecological criteria when selecting our global partners for storage and transportation. Each of our partners must offer a spectrum of services that clearly reduces greenhouse gases, from their fleets to their equipment.