

# Sartorius EHS Policy Statement

At Sartorius, we are committed to fostering safe, compliant workplaces and minimize the environmental impact of our business activities. The following principles guide our actions in EHS.



## Ensure Full Compliance

- Maintain an EHS Management System and implement robust processes to ensure all legal and internal standards are met while adapting quickly when necessary.
- Establish clear communication, and accountability.
- Ensure clear and comprehensive documentation and record keeping of all compliance-related actions and incidents.



## Prioritize Employee Health and Safety

- Never compromise on occupational health and safety for economic gain.
- Equip employees with comprehensive health and safety training.
- Actively develop programs to enhance physical, mental, and social well-being.
- Implementing rigorous safety measures and risk management to prevent accidents and health hazards.



## Drive EHS Excellence

- Foster a culture that aligns employee health and safety, environmental stewardship, and community well-being with our corporate values.
- Provide support for EHS goals, track progress, and conduct regular reviews.
- Share knowledge and best practices, and implement preventive actions to avoid incident reoccurrence.
- Encourage active participation in EHS groups, and report incidents and unsafe conditions promptly.



## Champion Environmental Protection

- Understand and act on our environmental impacts and uphold associated human rights.
- Provide targeted environmental training to our employees.



## Integrate EHS in Business

- Embed EHS into the early stages of product, process, and technological development, as well as procurement and investment decisions.



## Mobilize Business-Wide EHS Support

- Contribute to Sartorius' EHS and sustainability goals by working safely and responsibly.
- Strengthen EHS expertise through strategic hiring, training, and development initiatives.

Signiert von:

Joachim Kreuzburg

Name des Unterzeichners: Joachim Kreuzburg  
Signiergrund: Ich genehmige dieses Dokument  
Signierzeit: 2024-10-29 | 15:39:52 MEZ

EC96CF6C5CFB4391AC89FCB439227DA4

Signiert von:

Florian Presting

Name des Unterzeichners: Florian Presting  
Signiergrund: Ich habe dieses Dokument verfasst  
Signierzeit: 2024-10-29 | 15:04:32 MEZ

CAB21158101849F389259922421F6160

Dr. Joachim Kreuzburg  
CEO

Florian Presting  
Head of EHS & Security

Head of Site | Business