

FSC® Certified - Enhance Your Experience with Eco-Friendly Packaging Choices



What is FSC®?

The Forest Stewardship Council (FSC) is a global non-profit organization that promotes environmentally sound and socially beneficial management of the world's forests. FSC certification ensures that paper, packaging and wood-based materials are sourced sustainably from well-managed forests.

It certifies both forests and supply chains, guaranteeing compliance with high environmental and social standards. By choosing FSC-certified materials, companies support biodiversity preservation and sustainable resource use.

What FSC® Means for You?

The FSC label on a packaging or printed materials means:

- Reducing environmental impact
- Ensuring transparency across supply chain
- Supporting fair labor and community rights

Discover the Difference FSC® Makes



Environment & Climate

FSC® prevents deforestation and promotes responsible forest management.



United Nations Sustainable Development Goals

FSC® is an effective tool to help achieve 40 targets under the UN's 14 Sustainable Development Goals (SDGs).



Gender Equity & Workers' Rights

FSC® standards promote equality and workers' rights including gender equity.



Indigenous Peoples' Rights

FSC® protects Indigenous rights by requiring consent before forestry on traditional lands and ensuring their voices are heard.

At Sartorius, we are committed to delivering products in packaging that is both safe and easy for our customers to handle, while prioritizing sustainability. To achieve this, we annually evaluate the practices of our suppliers and the attributes of our packaging. This evaluation helps us track our progress towards zero deforestation goals and identify opportunities to increase the use of recycled and renewable materials. Additionally, we strive to enhance the percentage of packaging that carries sustainable forestry certifications, such as those awarded by the Forest Stewardship Council (FSC). The FSC® label assures that the materials used are sourced responsibly, with care for forests, respect for people, and consideration for future generations. We are proud to take another step toward a more sustainable future with the FSC® certification.

Discover how Sartorius supports your lab sustainability goals with innovative, eco-friendly solutions and FSC-certified packaging:

