

Goettingen, Germany,

Sartorius Statement on Conflict Minerals Regulation – June 2025

Dear Valued Customer,

On August 22, 2012, the U.S. Securities and Exchange Commission ("SEC") published regulations implementing Section 1502 of the Dodd-Frank Wall Street Reform and Consumer Protection Act ("Dodd-Frank Act"). The Dodd-Frank Act requires publicly traded companies to report their use of "Conflict Minerals" (Tantalum, Tin, Gold and Tungsten, or "3TG") to the U.S. Securities and Exchange Commission. While Sartorius AG and its subsidiaries ("Sartorius") are not US publicly traded companies, Sartorius is committed to complying with the Dodd-Frank Act and related laws and regulations which aim to prevent the use of Conflict Minerals that are directly or indirectly sourced from conflict-affected and high-risk areas.

Sartorius' commitment to conducting its business with respect for human rights and in compliance with all applicable laws is supported by Sartorius Values and Code of Conduct for Business Partners. Sartorius recognizes the risks of significant adverse impacts which may be associated with extracting, trading, handling and exporting minerals from conflict-affected and high-risk areas and Sartorius is committed to its social responsibility to respect human rights and not contribute to conflict. Based upon existing information, Sartorius does not knowingly use Conflict Minerals derived from the Conflict Region in its products.

Further, in support of Sartorius commitment, Sartorius expects its suppliers to continuously monitor both direct and indirect supply chains to avoid procurement of materials from conflict regions, and to be forthright in sharing compliance information with Sartorius. Sartorius suppliers are encouraged to exercise due diligence on the source and chain of custody of Conflict Minerals and shall be requested to make their due diligence measures available to Sartorius upon request.

For more information, please contact the responsible persons listed on Sartorius' Documents and Positions webpage found in www.sartorius.com.

Sincerely,

Signiert von:

Torsten Müller



Name des Unterzeichners: Torsten Müller
Signiergrund: Ich genehmige dieses Dokument
Signierzeit: 2025-06-11 | 15:58:07 MESZ
DB2E8BA700F247F1846FE7C91706C30B

Torsten Müller
Head of Operations BPS

Signiert von:

Carsten Wolter



Name des Unterzeichners: Carsten Wolter
Signiergrund: Ich genehmige dieses Dokument
Signierzeit: 2025-06-12 | 10:16:13 MESZ
714DF2C4E2EC4DEC9D70956252D2F540

Carsten Wolter
Head of Operations LPS