

Goettingen, Germany,

Sartorius Statement on Conflict Minerals Regulation – April 2023

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

DocuSigned by:
Eric Kummer
Nom du signataire : Eric Kummer
Motif de la signature : J'approuve ce document
Heure de signature : 2023-05-17 | 15:22:56 CEST
DCA82C960A124914A1C390D79AF86363

DocuSigned by:
Laurent Claisse
Signer Name: Laurent Claisse
Signing Reason: I approve this document
Signing Time: 2023-05-10 | 13:51:55 CEST
1EF8204DFB6344E98A058AAA8F9DECC2

Eric Kummer
Head of Operations BPS

Laurent Claisse
Head of Operations LPS