

November 22, 2024

Sartorius Tax Principles

Sartorius is committed to ensuring that the company's tax affairs are diligently managed and to be recognized by stakeholders as a responsible and reliable taxpayer. It is a matter of principle that Sartorius pays the taxes legally due in any country or territory, in accordance with the rules set by the respective governments. Sartorius recognizes that tax payments are an essential part of revenues for governments needed for the fulfillment of their social, economic and societal tasks. Sartorius complies with all relevant tax regulations as well as filing, documentation, disclosures and approval requirements.

1. Sartorius tax fundamentals

- Strong opposition against noncompliance and tax fraud
- Tax payments in line with respective local value creation
- Tax optimization in line with business needs; no engagement in artificial transactions without business substance and/or in profit-shifting activities
- If applicable and reasonable, usage of government-sponsored incentives and reliefs, such as tax credits
- Transparency and cooperative approach with tax authorities

2. Tax organization

To fulfill its tax duties, Sartorius has defined clear lines of accountability, responsibility and organization for the tax function: The regional Heads of Finance have responsibility for all taxes and tax effects relating to the company's activities within their regional responsibility and report regularly to the Head of Finance. Additionally, to ensure transactions at arm's length, all cross-border intercompany relationships are supervised by the Group Tax department which reports to the Head of Tax. The Head of Tax and the Head of Finance update the Chief Financial Officer on a regular basis on all important tax-related issues. Beyond, the Head of Finance reports regularly on tax matters to both the Executive Board and the Audit Committee of the Supervisory Board.

3. Tax compliance and risk management

Taxation of international business is complex, and uncertainties may arise in the application of tax laws, resulting in tax risks. Sartorius follows a balanced approach to tax risk, taking into account the economic and operational factors as well as social aspects and reputation. As part of our internal procedures the tax obligations are routinely monitored. Sartorius tax professionals are regularly trained and updated to ensure that the necessary knowledge and qualifications are available. Additionally, advice is also sought from third party advisors and tax specialists. Sartorius has set up a compliance management system and a risk reporting process driving tax compliance and limiting tax risks. Remaining transparent and having a cooperative relationship between Sartorius and the relevant tax authorities creates trust and reduces tax risks or disputes.

November 2024 | Approved by CEO/CFO