

## Sartorius Aktiengesellschaft Göttingen

## **Annual General Meeting 2024**

Information according to Table 8 of the Implementing Regulation (EU) 2018/1212

Type of information	Explanation
Unique identifier of the corporate event	SRT032024Div
2. Nature of the corporate event	Profit distribution
3. ISIN	DE0007165607
4. ISIN (other)	DE0007165631
5. Uniform Resource Locator (URL)	www.sartorius.com/agm
Further information	Distribution of a dividend of € 0.73 per ordinary share entitled to dividend for the past financial year 2023.
	Distribution of a dividend of € 0.74 per preference share entitled to dividend for the past financial year 2023.
	Ex-Day: 2024-04-02; Payday: as of 2024-04-04