

Sartorius Aktiengesellschaft
Göttingen

Annual General Meeting 2026

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

Type of information	Explanation
1. Unique identifier of the corporate event	SRT032026Div
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	<p>Distribution of a dividend of €0.73 per ordinary share entitled to dividend for the past financial year 2025.</p> <p>Distribution of a dividend of €0.74 per preference share entitled to dividend for the past financial year 2025.</p> <p>Ex-Day: 2026-03-27; Payday: as of 2026-03-31</p>