

Information Pursuant to Section 125(1) of the German Stock Corporation Act (AktG) in Conjunction with Section 125(5) of AktG and Article 4(1) as well as Table 3 of the Annex to Implementing Regulation (EU) 2018/1212

Type of Information	Description
A. Specification of the Message	
1. Unique identifier of the event	SRT032024oHV
2. Type of message	Invitation to the Annual General Meeting [in the format specified in Implementing Regulation (EU) 2018/1212: NEWM].
B. Specification of the Issuer	
1. ISIN	DE0007165607 and DE0007165631
2. Name of issuer	Sartorius Aktiengesellschaft
C. Specification of the Meeting	
1. Date of the General Meeting	March 28, 2024 [in the format specified in Implementing Regulation (EU) 2018/1212: 20240328].
2. Time of the General Meeting	12:00 p.m. (CET) [in the format specified in Implementing Regulation (EU) 2018/1212: 11:00 UTC]
3. Type of General Meeting	Annual General Meeting held as a virtual General Meeting without shareholders or their proxies being physically present [in the format specified in Implementing Regulation (EU) 2018/1212: GMET].
4. Location of the General Meeting	URL of the virtual General Meeting: www.sartorius.com/agm Location of the Annual General Meeting as required by the German Stock Corporation Act (AktG): Event hall at Annastr. 29, 37075 Göttingen, Germany
5. Record Date	The record date within the meaning of Sec. 123(4) of AktG is March 06, 2024 (midnight CET). [in the format specified in Implementing Regulation (EU) 2018/1212: 20240306, 23:00 UTC]
6. Uniform Resource Locator (URL)	www.sartorius.com/agm



D. Participation in the General Meeting D. Participation in the General Meeting – Voting by Electronic Communication		
	[in the format specified in Implementing Regulation (EU) 2018/1212: EV; ISO 20022: EVOT]	
2. Issuer deadline for the notification of participation	March 21, 2024 (midnight CET)	
	[in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC]	
3. Issuer deadline for voting	March 28, 2024, up to the time announced for this purpose by the chair of the meeting at the virtual Annual General Meeting	
	[in the format specified in Implementing Regulation (EU) 2018/1212:	
	20240328: up to the time announced for this purpose by the chair of the meeting at the virtual Annual General Meeting	
D. Participation in the General Meeting – Proxies Designated b	by the Company	
1. Method of participation by shareholder	Granting of power of attorney and instructions to the proxies designated by	
	the company [in the format specified in Implementing Regulation (EU) 2018/1212:	
	PX; ISO 20022: PRXY]	
2. Issuer deadline for the notification of participation	March 21, 2024 (midnight CET)	
	[in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC]	
3. Issuer deadline for voting	Granting of power of attorney (with instructions) to the proxies designated	
	by the company	
	 via postal mail or email by March 27, 2024, 6:00 p.m. (CET) [in the format specified in Implementing Regulation (EU) 2018/1212: 	
	20240327; 17:00 UTC]	
	 electronically via the online platform at www.sartorius.com/agm by March 28, 2024, by the time announced for 	
	this purpose at the Annual General Meeting by the chair of the meeting	
	[in the format specified in Implementing Regulation (EU) 2018/1212:	
	20240328; up to the time announced for this purpose by the chair of the meeting at the Annual General Meeting	
D. Participation in the General Meeting – Exercise of Voting Ri	ghts by an Authorized Third Party	
1. Method of participation by shareholder	Voting rights exercised by authorized third parties by means of electronic	
	communication or by granting sub-power of attorney and instructions to the proxies designated by the company	
	[in the format specified in Implementing Regulation (EU) 2018/1212:	
	PX; ISO 20022: PRXY]	
2. Issuer deadline for the notification of participation	March 21, 2024 (midnight CET)	
	[in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC]	
3. Issuer deadline for voting	Voting rights exercised by authorized third parties by means of electronic	
	communication via the online platform at www.sartorius.com/agm:	
	 March 28, 2024, , up to the time announced for this purpose by the chair of the meeting at the virtual Annual General Meeting 	
	[in the format specified in Implementing Regulation (EU) 2018/1212:	
	20240328: up to the time announced for this purpose by the chair of the meeting at the virtual Annual General Meeting	
	Voting rights exercised by authorized third parties by issuing sub-power of	
	attorney and instructions to the proxies designated by the company:	
	 via postal mail or email by March 27, 2024, 6:00 p.m. (CET) [in the format specified in Implementing Regulation (EU) 2018/1212: 	
	20240327; 17:00 UTC]	
	electronically via the online platform at www.sartorius.com/agm	
	by March 28, 2024, up to the time announced for this purpose at the Annual General Meeting by the chair of the meeting	
	Allitual Deficial Meetific by the Chall of the meetific	
	[in the format specified in Implementing Regulation (EU) 2018/1212: 20240328; up to the time announced for this purpose by the chair of the	



E. Agenda		
E. Agenda – Agenda Item 1		
1. Unique identifier of the agenda item	1	
2. Title of the agenda item	Presentation of the approved annual financial statements of Sartorius Aktiengesellschaft and the endorsed consolidated financial statements for the year ended December 31, 2023; the combined management report for Sartorius Aktiengesellschaft and the Group, together with the explanatory report, included therein, of the Executive Board concerning the disclosures according to Sections 289a and 315a of the German Commercial Code (HGB); as well as together with the report of the Supervisory Board for fisca 2023	
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm	
4. Vote		
5. Alternative voting options		
E. Agenda - Agenda Item 2		
1. Unique identifier of the agenda item	2	
2. Title of the agenda item	Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft	
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm	
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]	
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]	
E. Agenda - Agenda Item 3		
1. Unique identifier of the agenda item	3	
2. Title of the agenda item	Resolution on granting discharge to the members of the Executive Board for fiscal 2022	
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm	
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]	
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]	
E. Agenda - Agenda Item 4		
1. Unique identifier of the agenda item	4	
2. Title of the agenda item	Resolution on granting discharge to the members of the Supervisory Board for fiscal 2023	
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm	
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]	
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]	



E. Agenda - Agenda Item 5	<u>_</u>
Unique identifier of the agenda item	5
2. Title of the agenda item	Appointment of an auditor for fiscal 2024 as well as an auditor for the audit review or audit of interim financial reports financial information in fiscal 2024 and in fiscal 2025 in the period up to the next Annual General Meeting
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]
E. Agenda – Agenda Item 6	
1. Unique identifier of the agenda item	6
2. Title of the agenda item	Resolution on approval of the remuneration report
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm
4. Vote	Advisory vote [in the format specified in Implementing Regulation (EU) 2018/1212: AV; ISO 20022: ADVI]
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]
E. Agenda – Agenda Item 7	
1. Unique identifier of the agenda item	7
2. Title of the agenda item	Resolution on granting a new authorization pursuant to Section $71(1)(8)$ of the German Stock Corporation Act (AktG) relating to the acquisition and the use of treasury shares with the option of excluding subscription rights
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]
E. Agenda - Agenda Item 8	
1. Unique identifier of the agenda item	8
2. Title of the agenda item	Resolution on an authorization to use derivatives in connection with the acquisition of treasury shares with the exclusion of shareholders' subscription and tender rights
3. Uniform Resource Locator (URL) of the materials	www.sartorius.com/agm
4. Vote	Binding vote [in the format specified in Implementing Regulation (EU) 2018/1212: BV; ISO 20022: BNDG]
5. Alternative voting options	Vote in favor, vote against, abstention [in the format specified in Implementing Regulation (EU) 2018/1212: VF, VA, AB; ISO 20022: CFOR, CAGS, ABST]



F. Shareholders' Rights – Additions to the Agenda Pursuant to Section 122(2) of AktG		
1. Object of deadline	Submission of the request to add items to the agenda	
2. Applicable issuer deadline	February 26, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240226; 23:00 UTC]	
F. Shareholders' Rights – Countermotions Pursuant	to Section 126(1) and (4) of AktG	
1. Object of deadline	Submission of countermotions to specific agenda items	
2. Applicable issuer deadline	March 13, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240313; 23:00 UTC]	
F. Shareholders' Rights – Election Proposals Pursuar	nt to Section 127 of AktG	
1. Object of deadline	Submission of proposals for the election of Supervisory Board members or auditors	
2. Applicable issuer deadline	March 13, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240313; 23:00 UTC	
F. Shareholders' Rights – Submission of Statements	Pursuant to Section 130a(1) to (4) of AktG	
1. Object of deadline	Submission of statements on agenda items by means of electronic communication	
2. Applicable issuer deadline	For registering for the Annual General Meeting: March 21, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC] For submitting statements: March 22, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240322; 23:00 UTC]	
	ing Pursuant to Section 130a(5) of AktG (including the right to submit motions and election and all types of requests for information pursuant to Section 131 of AktG, in each case as part	
1. Object of deadline	Right to speak at the meeting by way of video communication. Motions and election proposals in accordance with Section 118a(1)(2)(3) of AktG as well as all types of requests for information in accordance with Section 131 of AktG may be part of the address. It is intended that the chair of the virtual Annual General Meeting will stipulate in accordance with section 131(1f) of AktG that the right to information pursuant to section 131 of AktG may only be exercised by mear of video communication.	
2. Applicable issuer deadline	For participating in the Annual General Meeting: March 21, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC] For exercising the right to speak: March 28, 2024; during the Annual General Meeting as specified by the chair of the meeting [in the format specified in Implementing Regulation (EU) 2018/1212: 20240328; during the Annual General Meeting as specified by the chair of the meeting]	



F. Shareholders' Rights – Submission of the Request Pursuant to Section 131(4)(1) and (2) of AktG		
1. Object of deadline	Submission of the request pursuant to Section 131(4)(1) and (2) of AktG by means of electronic communication	
2. Applicable issuer deadline	For participating in the Annual General Meeting: March 21, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC] For transmitting the request: March 28, 2024; during the Annual General Meeting until the time determined by the chair of the meeting [in the format specified in Implementing Regulation (EU) 2018/1212: 20240328; during the Annual General Meeting until the time determined by the chair of the meeting]	
F. Shareholders' Rights – Submission of the Reques	t Pursuant to Section 131(5)(1) and (2) of AktG	
1. Object of deadline	Submission of the request pursuant to Section 131(5)(1) and (2) of AktG by means of electronic communication	
2. Applicable issuer deadline	For participating in the Annual General Meeting: March 21, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC] For transmitting the request: March 28, 2024; until the Annual General Meeting is closed by the chair of the meeting [in the format specified in Implementing Regulation (EU) 2018/1212: 20240328; until the Annual General Meeting is closed by the chair of the meeting]	
F. Shareholders' Rights - Objection		
1. Object of deadline	Submitting objections to resolutions of the Annual General Meeting by means of electronic communication	
2. Applicable issuer deadline	For participating in the Annual General Meeting: March 21, 2024, midnight (CET) (time of receipt by the company) [in the format specified in Implementing Regulation (EU) 2018/1212: 20240321; 23:00 UTC] For submitting objections: on March 28, 2024 from the beginning of the virtual Annual General Meeting until the Annual General Meeting is closed by the chair of the meeting. [in the format specified in Implementing Regulation (EU) 2018/1212: 20240328; from the beginning of the virtual Annual General Meeting until the Annual General Meeting is closed by the chair of the meeting]	

Other Disclosures Pursuant to Section 125(1)(4) and (5) of AktG

Voting rights may be exercised by proxy, including by a union of shareholders

. . .

Further disclosures can also particularly be found in the invitation to the Annual General Meeting, which is available on the company's website at www.sartorius.com/agm.