

Annual General Meeting of Sartorius Aktiengesellschaft on March 26, 2026

Granting of Power of Attorney and Issuing Instructions to Proxies Designated by the Company

ID number confirming registration (AGM ticket): _____ Last name | Company: _____
Number of ordinary shares: _____ First name: _____
Phone number*: _____ Email address*: _____

Please return by no later than March 25, 2026, 6:00 p.m. (CET) (deadline for receipt by the company) to:

Sartorius Aktiengesellschaft or by **email**: sartorius@captrace.com
c/o CAPTRACE GmbH
Trimburgstr. 2
81249 Munich
Germany

I | we grant power of attorney for the virtual general meeting on March 26, 2026, to each of the proxies designated by the company, Ms. Carolin Brandt and Mr. Jens M. Artelt, both employees of Sartorius Corporate Administration GmbH that is legally domiciled in Göttingen, Germany, and who are based in Göttingen, Germany, each individually and with the right to issue substitute powers of attorney, to represent me | us and to exercise my | our voting right(s) as marked by an X on the next page. Power of attorney and instructions previously issued to the proxies designated by the company, as well as voting by (electronic) communication, are hereby revoked.

I | we hereby confirm that I | we have read the information on granting a power of attorney and instructions on voting to the proxies designated by the company and accept the conditions on the proxy voting listed therein.

 Be sure to send this page and the next one.

Please go to the next page to issue instructions on voting on the items of the agenda

Instructions on Proxy Voting

The instructions each relate to the proposed resolutions from the company's boards for the individual items on the agenda announced on February 16, 2026, in the German Federal Gazette (Bundesanzeiger), if and to the extent that no instructions are given relating to counter motions and | or nominations from shareholders (in relation to this, please refer to the explanations at the end of this document). The instructions relating to the resolution on agenda item 2 (resolution on the appropriation of the retained profit) proposed by the company's boards likewise apply to a proposed resolution by the company's boards that is put to the vote at the Annual General Meeting in a correspondingly adjusted form, if necessary, as a result of a change in the number of shares carrying dividend rights, provided that the instructions or votes are not amended or revoked. If an individual vote is held on an agenda item instead of a collective vote without this having been communicated in advance of the Annual General Meeting, the instructions issued to the proxies on this agenda item will apply accordingly to each item of the associated individual vote, unless they are amended or revoked. Please mark the "Yes" field with an X if you agree, or the "No" field with an X if you reject the resolution. If no vote should be cast (abstention), please mark the "Abstain" field with an X.

Agenda item:

	Yes	No	Abstain
2. Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
3. Resolution on granting discharge to the members of the Executive Board for fiscal year 2025	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
4. Resolution on granting discharge to the members of the Supervisory Board for fiscal year 2025	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
5.1 Appointment of the auditor for fiscal year 2026 as well as the auditor for the audit review or audit of interim financial reports financial information in fiscal year 2026 and in fiscal year 2027 in the period up to the next Annual General Meeting	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
5.2 Appointment of the auditor for the sustainability report for fiscal year 2026	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
6. Resolution on approval of the remuneration report	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>

Counter motions and Nominations from Shareholders

Any counter motions and nominations from shareholders received by midnight (CET) on March 11, 2026, will be published online at www.sartorius.com/aggm. Information can also be found there about how you can join counter motions and nominations. If you wish to issue instructions on proxy voting for the relevant counter motions | nominations, please place an X in the "Yes" field if you agree, or the "No" field if you disagree. If you do not wish to cast a vote (abstention), please place an X in the "Abstain" field.

Motion Nomination	Yes	No	Abstain	Motion Nomination	Yes	No	Abstain
A	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	D	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
B	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	E	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
C	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	F	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>

Place

Date

Legible signature(s) or person(s) making this declaration

Instructions on Granting Power of Attorney and Issuing Instructions to the Proxies Designated by Sartorius Aktiengesellschaft

You may grant power of attorney to the proxies who are designated by Sartorius Aktiengesellschaft to vote on your behalf. These proxies, Ms. Carolin Brandt and Mr. Jens M. Artelt, both employees of Sartorius Corporate Administration GmbH that is legally domiciled in Göttingen, Germany, and who are based in Göttingen, Germany, are bound by your voting instruction, and each shall have sole power of representation. As a result of your power of attorney, the proxies will be entitled to vote on your behalf only if you have given them **explicit instructions** on how they are to exercise the voting rights. The proxies are obligated to vote according to your instructions. You have the following options for granting power of attorney and issuing instructions to the proxies mentioned above using the form “Granting Power of Attorney and Issuing Instructions to Proxies Designated by the Company”:

Sending of the completed form for granting power of attorney and issuing instructions to the proxies designated by Sartorius Aktiengesellschaft by postal mail or email

For this purpose, please use the form “Granting Power of Attorney and Issuing Instructions to Proxies Designated by the Company.” Grant power of attorney to the aforementioned proxies of Sartorius Aktiengesellschaft and instruct them how to exercise your voting rights on the resolutions proposed by the company’s boards by marking the appropriate boxes with an X. Then send your “Granting Power of Attorney and Issuing Instructions to Proxies Designated by the Company” form by postal mail or email, indicating your AGM ticket number confirming your registration, directly to the following address by no later than March 25, 2026, 6:00 p.m. (CET) (deadline for receipt by the company):

- By postal mail to:
Sartorius Aktiengesellschaft
c/o CAPTRACE GmbH
Trimbургstr. 2
81249 Munich
Germany
- By email to: sartorius@captrace.com

Alternatively, you can use the AGM Portal available on the company’s website at www.sartorius.com/agm by no later than the time announced in the virtual Annual General Meeting by the chair of the meeting for the respective vote to grant power of attorney and issue instructions to proxies designated by the company in accordance with the procedure defined there.

Important information:

Please note that only after registering in due form and time are shareholders entitled to grant powers of attorney and issue voting instructions to the proxies designated by Sartorius Aktiengesellschaft. You can send your completed form along with your confirmation of registration or submit your AGM ticket ID number no later than March 25, 2026, 6:00 p.m. (CET) (deadline for receipt by the company). If the company receives divergent declarations on the exercise of voting rights for the same shareholding by different means of transmission, only the declaration received last will be taken into account. The appointment of proxies designated by the company to exercise other shareholder rights, in particular to submit countermotions, ask questions and to file objections, is excluded. Even after you have granted power of attorney to the proxies designated by the company, you are still entitled to use the AGM Portal with all the options it offers, according to the procedure it provides for this purpose, up until the end of the virtual general meeting on March 26, 2026.

If you have any questions regarding voting by proxy, please contact our **Annual General Meeting hotline** from Monday to Friday, except public holidays, between 9:00 a.m. and 5:00 p.m. at **+49 (0)611 711 868 49**.