

Annual General Meeting of Sartorius Aktiengesellschaft on March 29, 2023

Granting of Power of Attorney and Issuing Instructions to Voting 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 28, 2023, 6:00 p.m. (CEST)** (deadline for receipt by the company) to:

Sartorius Aktiengesellschaft
c/o Better Orange IR & HV AG
Haidelweg 48
81241 Munich
Germany

or by **email**: sartorius@better-orange.de

I | we grant power of attorney for the virtual general meeting on March 29, 2023, to each of the voting proxies designated by the company, Ms. Katrin Sebastian and Mr. Jens M. Artelt, both authorized signatories of Sartorius Aktiengesellschaft 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 below. Power of attorney and instructions previously issued to the voting proxies designated by the company, as well as voting by (electronic) correspondence, 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 voting 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 management for the individual items on the agenda announced on February 17, 2023, in the German Federal Gazette (Bundesanzeiger), if and to the extent that no instructions are given relating to countermotions and | or nominations from shareholders (in relation to this, please refer to the explanations at the end of this page). The instructions relating to the resolution on agenda item 2 (resolution on the appropriation of retained profits) proposed by the company's management bodies likewise apply to a proposed resolution by the company's management bodies 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, insofar as the instructions 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 2022	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
4. Resolution on granting discharge to the members of the Supervisory Board for fiscal 2022	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
5. Appointment of an auditor for fiscal 2023 as well as an auditor for the audit review of the first-half financial report of 2023	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
6. Resolution on the approval of the remuneration policy for the Executive Board members	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
7. Resolution on approval of the remuneration report	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
8. Resolution on an amendment to the remuneration paid to the members of the Supervisory Board, a corresponding amendment to Article 12 of the Articles of Association (Remuneration) and the associated system for the remuneration of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
9. Resolution on an amendment to the Articles of Association to enable virtual annual general meetings by adding a new Article 13a to the Articles of Association	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
10. Resolution on an amendment to Article 15 of the Articles of Association (Entitlement to Attend) concerning the participation of Supervisory Board members in the Annual General Meeting by means of video and audio communication	<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 14, 2023, will be published online at www.sartorius.com/agm. 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) of person(s) making this declaration

Instructions on Granting Power of Attorney and Issuing Instructions to the Voting 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 voting proxies, Ms. Katrin Sebastian and Mr. Jens M. Artelt, both authorized signatories of Sartorius Aktiengesellschaft that is legally domiciled in Göttingen, Germany, and who are based in Göttingen, Germany, are bound by your voting instructions; and each shall have sole power of representation. As a result of your power of attorney, the voting proxies will be entitled to vote on your behalf only if you have given them **explicit instructions** on how they are to vote on the individual resolution proposals of company management in the agenda items.

The voting proxies are obligated to vote according to your instructions on the resolution proposals of company management announced in the agenda. You have the following options for granting power of attorney and issuing instructions to the voting proxies mentioned above using the form "Granting Power of Attorney and Issuing Instructions to Voting Proxies Designated by the Company":

Sending of the completed form for granting power of attorney and issuing instructions to the voting 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 Voting 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 management bodies. Then send your "Granting Power of Attorney and Issuing Instructions to Voting 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 28, 2023, at 6:00 p.m. (CEST) (deadline for receipt by the company):

- By mail (post) addressed to: Sartorius Aktiengesellschaft
c/o Better Orange IR & HV AG
Haidelweg 48
81241 Munich
Germany
- via email to: sartorius@better-orange.de

Alternatively, you can use the password-protected online platform for the Annual General Meeting available on the company's website at www.sartorius.com/agm up to the time announced by the chair of the virtual Annual General Meeting to grant power of attorney and issue instructions to voting 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 voting 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 28, 2023, at 6:00 p.m. (CEST) (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 voting proxies designated by the company to submit countermotions, ask questions and to file objections is excluded. Even after you have granted power of attorney to the voting proxies designated by the company, you are still entitled to use the password-protected Internet service platform 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 29, 2023.

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 **+49-89-889-690-6610**.