

Annual General Meeting of Sartorius Aktiengesellschaft on March 28, 2024

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 adress*:
Please return by no later than March 27, 2024, 6:00 p.	m. (CET) (deadline for receipt by the company) to:
Sartorius Aktiengesellschaft or by email : sa c/o Better Orange IR & HV AG Haidelweg 48 81241 Munich Germany	artorius@better-orange.de

I | we grant power of attorney for the virtual general meeting on March 28, 2024, 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

*Voluntary information 1



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 16, 2024, 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 document). The instructions relating to the resolution on agenda item 2 (resolution on the appropriation of retained profits) proposed by the company's management likewise apply to a proposed resolution by the company's management 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	N	10	Abstain	
2.	. Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft									
3.	 Resolution on granting discharge to the members of the Executive Board for fiscal 2023 									
4.										
5.	5. 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 MeetingResolution on approval of the remuneration report							Г	\neg		
7.								=		
	of the German Stock Corporation Act (AktG) relating to the acquisition and the							_		
Q	use of treasury shares with the option of excluding subscription rights Resolution on an authorization to use derivatives in connection with the acquisition						Г	\neg		
8. Resolution on an authorization to use derivatives in connection with the acquisition of treasury shares with the exclusion of shareholders' subscription and tender rights										
Countermotions and Nominations from Shareholders Any countermotions and nominations from shareholders received by midnight (CET) on March 13, 2024, will be published online at www.sartorius.com/agm. Information can also be found there about how you can join countermotions and nominations. If you wish to issue instructions on proxy voting for the relevant countermotions 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.										
Мо	tion Nomination	Yes	No	Abstain	Motion Nomination	Yes	No	Abstain	1	
Α					D					
В					Е					
С					F					
——Plac	e		te.		egible signature(s) of person(s)	making t	his declar	ation		



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 management 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 27, 2024, 6:00 p.m. (CET) (deadline for receipt by the company):

■ By postal mail to: Sartorius Aktiengesellschaft

c/o Better Orange IR & HV AG

Haidelweg 48 81241 Munich Germany

By email to: sartorius@better-orange.de

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 27, 2024, 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 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 28, 2024.

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)89 889 690 6610**.