

# Virtual General Meeting of Sartorius Aktiengesellschaft on June 26, 2020

Voting by Mail (Postal Voting) | 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 **no later than June 25, 2020, at 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  
or by **fax**: +49 (0)89 889 690 655

## Please clearly mark the appropriate boxes with an X:

(If an X is missing or a box is not clearly marked, voting by mail shall be assumed. If several declarations of intent are received, the point in time at which these are received shall determine which version shall count. If the company receives declarations differing from one another for one and the same share through different transmission channels and if it is not possible for the company to determine which declaration was last submitted, these declarations will be treated as binding in the following order of transmission channels: (1) Internet service platform for the virtual general meeting, (2) email, (3) fax and (4) paper form.)

- I | We exercise my | our voting rights with respect to the virtual general meeting on June 26, 2020, by postal vote as marked below. Votes cast earlier are hereby revoked.
- I | we grant power of attorney 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, to represent me | us and to exercise my | our voting right(s) as marked by an X below during the virtual general meeting on June 26, 2020. Power of attorney and instructions previously issued to the voting proxies designated by the company are hereby revoked.

Please go to the next page to cast votes by way of voting by mail or to issue instructions on voting on the items of the agenda

Cast votes by way of voting by mail on the items of the agenda or issue voting instructions for these

	Yes <sup>1</sup>	No
2. Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft	<input type="checkbox"/>	<input type="checkbox"/>
3. Resolution on granting discharge to the members of the Executive Board for fiscal 2019	<input type="checkbox"/>	<input type="checkbox"/>
4. Resolution on granting discharge to the members of the Supervisory Board for fiscal 2019	<input type="checkbox"/>	<input type="checkbox"/>
5. Resolution on the amendment of the Articles of Association	<input type="checkbox"/>	<input type="checkbox"/>
6. Elections for the Supervisory Board: elect Prof. David Ebsworth	<input type="checkbox"/>	<input type="checkbox"/>
7. Appointment of an auditor for fiscal 2020 as well as an auditor for the audit review of the first-half financial report of 2020	<input type="checkbox"/>	<input type="checkbox"/>

\_\_\_\_\_

Place

\_\_\_\_\_

Date

\_\_\_\_\_

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

**Please be sure to send both completed pages!**

## Instructions on Voting by Mail | Granting Power of Attorney and Issuing Instructions to the Voting Proxies Designated by Sartorius Aktiengesellschaft

You may vote by mail (post) or 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, are both authorized signatories of Sartorius Aktiengesellschaft that is legally domiciled in Göttingen, Germany; are based in Göttingen, Germany; will be bound by your voting instructions; and each shall have sole power of representation. 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 (“the company administration”) regarding the agenda. The voting proxies are obligated to vote according to your instructions on the resolution proposals announced in the agenda. You have the following options for voting by mail and for granting power of attorney to the voting proxies mentioned above using the form “Voting by Mail (Postal Voting) | Granting Power of Attorney and Issuing Instructions to Voting Proxies Designated by the Company”:

### **Sending by post, faxing or emailing the completed form for voting by mail or for granting power of attorney and issuing instructions to the voting proxies designated by Sartorius Aktiengesellschaft**

For this purpose, please use the form “Voting by Mail (Postal Voting) | Granting Power of Attorney and Issuing Instructions to Voting Proxies Designated by the Company.” Vote by mail or grant power of attorney to the above-mentioned proxies designated by Sartorius Aktiengesellschaft and issue instructions to these persons on how to vote on the resolution proposals of company management (“the company administration”). Then send this form by post, fax or email, indicating your AGM ticket number confirming your registration, directly to the following address:

- By mail (post) addressed to: Sartorius Aktiengesellschaft  
c/o Better Orange IR & HV AG  
Haidelweg 48  
81241 Munich, Germany
- Or by fax sent to the following number: +49 (89) 889 690 655
- Or by email sent to: sartorius@better-orange.de

**As an alternative, you can use the password-protected Internet service platform for the Annual General Shareholders’ Meeting (virtual general meeting) on the company’s Internet site at [www.sartorius.com/shareholders-meeting](http://www.sartorius.com/shareholders-meeting) according to the procedure provided there.**

### **Important information:**

Please note that only after registering in due form and time are shareholders entitled to exercise their voting rights by mail and 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 June 25, 2020, at 6:00 p.m. (CEST) (deadline of receipt by the company). If the company receives both a form for voting by mail and a power of attorney and instructions for the voting proxies designated by the company for one and the same ordinary share, the completed form for voting by mail shall solely be considered in any case. Furthermore, if the company receives declarations that differ from one another in connection with the granting and revocation of a power of attorney or proxy instructions via different transmission channels, and if it is not possible for the company to determine which of these declarations was last made, these declarations will be treated as binding in the following order of transmission: (1) Internet service platform for the Annual General Shareholders’ Meeting (virtual general meeting), (2) email, (3) fax and (4) paper form. The appointment of voting proxies designated by the company to submit countermotions, ask questions and to file objections is excluded. Even after you have voted by mail or 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 June 26, 2020.

If you have any questions about voting by mail (postal voting) and granting power of attorney to proxies, the staff members of our **Annual General Shareholders’ Meeting hot line** will be available to help you on Mondays through Fridays, except for legal holidays, between 9:00 a.m. and 5:00 p.m. (CEST) at **+49 (89) 889 690 620**.