

Virtual General Meeting of Sartorius Aktiengesellschaft on June 26, 2020

## Revocation of Voting by Mail

## Revocation of Power of Attorney Granted to a Third Party

## Revocation of Power of Attorney Granted and Instructions Issued to the Voting Proxies Designated by the Company

ID number confirming registration (AGM ticket): \_\_\_\_\_ Last name | Company: \_\_\_\_\_

Number of 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 box(es) with an X:**

I | we hereby **revoke** my | our votes cast **by mail** for the virtual general meeting on June 26, 2020.

I | we hereby **revoke power of attorney** granted **to the voting proxies designated by the company** to represent me | us during the virtual general meeting on June 26, 2020.

I | we hereby **revoke power of attorney** granted to a third party named below to represent me | us during the virtual general meeting on June 26, 2020.

Last name: \_\_\_\_\_

First name: \_\_\_\_\_

Place of residence: \_\_\_\_\_

Furthermore, I | we draw the former proxy's or company-appointed proxies' attention to the special condition that he | she is or they are **not permitted to use** the individual access data I | we have disclosed to him/her for the password-protected Internet service platform.

\_\_\_\_\_  
Place Date Legible signature(s) of person(s) making this declaration

**Please note the following:**

- Revocation of power of attorney must be in the form of text (§ 126b of the German Civil Code, a legible declaration on a tamper-proof permanent data medium that is inalterable once recorded) or must be submitted using the company's input form provided on the Internet service platform for the Annual General Shareholders' Meeting (virtual general meeting) at **[www.sartorius.com/shareholders-meeting](http://www.sartorius.com/shareholders-meeting)**.
- If you would like to revoke power of attorney granted to the voting proxies designated by the company, please send this revocation by June 25, 2020, 6:00 p.m. (CEST), to the following address:
- Sartorius Aktiengesellschaft, c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munich, Germany; or by fax: +49 (89) 889 690 655; or by email to **[sartorius@better-orange.de](mailto:sartorius@better-orange.de)**.
- After June 25, 2020, 6:00 p.m. (CEST), you will be able to revoke this power of attorney only by using the password protected Internet service platform for the virtual general meeting on the company's website at **[www.sartorius.com/shareholders-meeting](http://www.sartorius.com/shareholders-meeting)** according to the procedure provided there for this purpose.
- If you would like to revoke power of attorney granted to a third party, you may declare this revocation either to the third party you appointed as your proxy or to the company. In the event you wish to declare revocation of this power of attorney granted to a third party with respect to the company, please send this revocation by June 25, 2020, 6:00 p.m. (CEST), to the following address: Sartorius Aktiengesellschaft, c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munch, Germany; or by fax: +49 (89) 889 690 655; or by email to **[sartorius@better-orange.de](mailto:sartorius@better-orange.de)**.
- After June 25, 2020, 6:00 p.m. (CEST), you will be able to declare revocation of this power of attorney to the company only by using the password protected Internet service platform for the virtual general meeting on the company's website at **[www.sartorius.com/shareholders-meeting](http://www.sartorius.com/shareholders-meeting)** according to the procedure provided there for this purpose.