Annual General Meeting of Sartorius Aktiengesellschaft on March 28, 2024

Revocation of Power of Attorney Granted to a Third Party

Revocation of Power of Attorney Granted and Instructions Issued to the 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 by no later than March 27, 2024, 6:00 p.m. (CET) (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

Please clearly mark the appropriate box with an X:

	oke power of attorney granted to the proxies designated by al General Meeting on March 28, 2024.	<b>y the company</b> to represent me   us
	<b>yoke power of attorney granted to the third party named b</b> Meeting on March 28, 2024.	<b>below</b> to represent me   us during the
Last name:		
First name:		
Place of residence:		

Furthermore, I | we draw the former proxy's attention to the special condition that he | she **is not permitted to use** the individual access data that he | she has received for the AGM Portal.

## SVIFCTFAS

## Please note the following important information:

- Revocation of power of attorney must be submitted in text form (section 126b of the German Civil Code (BGB)) or, alternatively, is possible using the form on the company's AGM Portal available 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.
- If you would like to revoke the power of attorney granted to the proxies designated by the company, please send this revocation by March 27, 2024, 6:00 p.m. (CET), to the following address: Sartorius Aktiengesellschaft, c/o Better Orange IR & HVAG, Haidelweg 48, 81241 Munich, Germany, or via email to sartorius@better-orange.de.
- After March 27, 2024, 6:00 p.m. (CET), you will only be able to revoke this power of attorney via the AGM Portal available on the company's website at www.sartorius.com/agm in accordance with the procedure defined there.
- If you would like to revoke a 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 that you would like to revoke this power of attorney granted to a third-party proxy by notifying the company, please send this revocation to the following address by March 27, 2024, 6:00 p.m. (CET): Sartorius Aktiengesellschaft, c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munich, Germany, or via email to sartorius@better-orange.de.
- After March 27, 2024, 6:00 p.m. (CET), you will only be able to notify the company of your revocation of this power of attorney via the AGM Portal available on the company's website at www.sartorius.com/agm in accordance with the procedure defined there.