

Annual General Meeting of Sartorius Aktiengesellschaft on March 28, 2024

## Power of Attorney Granted to a Third Party

ID number confirming registration (AGM ticket):		Last name   Company:
registration (AON ticket).		Company.
Number of shares:		First name:
Phone number*:		Email address*:
Please return by <b>no later thar</b>	n March 27, 2024, 6:00 p	o.m. (CET) (deadline for receipt by the company) to:
Sartorius Aktiengesellschaft c/o Better Orange IR & HV AC Haidelweg 48 81241 Munich, Germany	_	sartorius@better-orange.de
I   we hereby grant power of a	attorney to Mr.   Mrs.	
Last name of proxy:		First name of proxy:
Proxy's place of residence:		
communication, to represent Meeting of Sartorius Aktiengename(s). This power of attorn	t me   us in exercising my esellschaft on March 28, ey includes the right of g	granted at an earlier time and any earlier voting by (electronic) our shareholder rights with respect to the virtual Annual General 2024, and to exercise my our voting rights by disclosing my our ranting substitute power of attorney. I we point out to said proxy she may not pass on said data to third parties.
Place	Date	Legible signature(s) of person(s) making this declaration

**Important information:** After receipt of power of attorney or proof of power of attorney granted, individual access data for the AGM Portal will be generated for the proxy. If the address details of the proxy are complete, the access data will be sent directly to the proxy along with a separate AGM ticket. If the address details of the proxy are incomplete, the AGM ticket for the proxy along with the individual access data for the AGM Portal will be sent to the grantor of the power of attorney for forwarding to the proxy. Please allow for usual processing and postal delivery times for the transmission of the access data. Special requirements may apply for granting power of attorney to banks or other intermediaries, unions of shareholders, proxy advisors or to other persons or institutions equal to the same pursuant to Section 135(8) of the German Stock Corporation Act (AktG). For this reason, shareholders are asked to consult with the proxies concerned, especially concerning the particular form of granting them power of attorney.

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 by the chair of the virtual Annual General Meeting for the respective vote to grant power of attorney to a third party in accordance with the procedure defined there.