

Annual General Meeting of Sartorius Aktiengesellschaft on March 26, 2026

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 25, 2026, 6:00 p.m. (CET)** (deadline for receipt by the company) to:

Sartorius Aktiengesellschaft or by **email**: sartorius@captrace.com
c/o CAPTRACE GmbH
Trimbургstr. 2
81249 Munich, Germany

Please clearly mark the appropriate box with an X:

I | we hereby revoke power of attorney granted to the proxies designated by the company to represent me | us during the Annual General Meeting on March 26, 2026.

I | we hereby revoke power of attorney granted to the third party named below to represent me | us during the Annual General Meeting on March 26, 2026.

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.

_____ Place

_____ Date

_____ Legible signature(s) or person(s) making this declaration

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**.
- If you would like to revoke the power of attorney granted to the proxies designated by the company, please send this revocation by March 25, 2026, 6:00 p.m. (CET), to the following address: Sartorius Aktiengesellschaft, c/o CAPTRACE GmbH, Trimbургstr. 2, 81249 Munich, Germany, or via email to **sartorius@captrace.com**.
- After March 25, 2026, 6:00 p.m. (CET), you will only be able to revoke the 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 a power of attorney granted to a third-party proxy by notifying the company, please send this revocation to the following address by
- March 25, 2026, 6:00 p.m. (CET): Sartorius Aktiengesellschaft, c/o CAPTRACE GmbH, Trimbургstr. 2, 81249 Munich, Germany, or via email to **sartorius@captrace.com**.
- After March 25, 2026, 6:00 p.m. (CET), you will only be able to notify the company of your revocation of the 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.

For further details, please refer to the information in the invitation to the Annual General Meeting.