

Annual General Meeting of Sartorius Aktiengesellschaft on March 26, 2026

Power of Attorney Granted to a Third Party

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

I | we hereby grant power of attorney to Mr. | Mrs.

Last name of proxy: _____ First name of proxy: _____

Proxy's place of residence: _____

if necessary by revoking any power of attorney already granted at an earlier time and any earlier voting by (electronic) communication, to represent me | us in exercising my | our shareholder rights with respect to the virtual Annual General Meeting of Sartorius Aktiengesellschaft on March 26, 2026, and to exercise my | our voting rights by disclosing my | our name(s). This power of attorney includes the right of granting substitute power of attorney. I | We point out to said proxy that upon receiving his | her individual access data to the AGM Portal, he | she may not pass on said data to third parties.

Place	Date	Legible signature(s) or 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 consider 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, associations 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 regarding the details, in particular the respective form of the power of attorney.

Alternatively, you can use the AGM Portal available on the company's website at www.sartorius.com/agm up to the end of the Annual General Meeting to grant power of attorney to a third party in accordance with the procedure defined there. If the proxy is to exercise the shareholder's voting rights, it is necessary to ensure that power of attorney is granted in good time before the vote in question.