

Annual General Shareholders' Meeting on March 28, 2019



Revocation of Power of Attorney

for entrance card no. _____ for _____ no-par shares in Sartorius Aktiengesellschaft

I/we _____ hereby revoke the power of attorney I/we granted

on _____ 2019 to

the voting proxies designated by Sartorius Aktiengesellschaft, Ms. Sabrina Hille and Mr. Mirko Koch, both residents of Göttingen, Germany,

Mr./Ms./Mrs. _____
Authorized third party

resident in _____
City

to represent me/us at the Annual General Shareholders' Meeting of Sartorius Aktiengesellschaft convened for **March 28, 2019**, and to exercise my/our rights (in particular voting rights).

Place/date Signature

Notes:

Revocations of powers of attorney must be made in the form of text (§ 126b of the German Civil Code – BGB).

Should you wish to revoke a power of attorney granted to the voting proxies designated by the company, please forward your revocation no later than by the end of **March 27, 2019 (midnight, 24:00 hours Central European Winter Time (CEWT), time of receipt)** to the following address:

Sartorius Aktiengesellschaft c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munich, Germany,
Fax: +49(0)89.889.690.633, Email: sartorius@better-orange.de

After this time, we would request that you hand in your revocation of the power of attorney at the entrance desk at the Annual General Shareholders' Meeting.

Should you wish to revoke power of attorney granted to a third party, you may notify said third party authorized of such revocation or alternatively notify the company as well. Should you wish to notify the company of such revocation, please forward your revocation no later than by the end of **March 27, 2019 (midnight, 24:00 hours CEWT, time of receipt)** to the following address:

Sartorius Aktiengesellschaft c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munich, Germany,
Fax: +49(0)89.889.690.633, Email: sartorius@better-orange.de

After this time, we would request that you hand in your revocation of the power of attorney at the entrance desk at the Annual General Shareholders' Meeting.

Sartorius Aktiengesellschaft