

Annual General Shareholders' Meeting on April 7, 2016



Revocation of Power of Attorney

for entrance card no. _____ for _____ ordinary shares in Sartorius AG

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

on _____ 2016 to

☐ the voting proxies designated by Sartorius AG, Ms. Inga Stucke and
Ms. Marina Ciottariello, both residents of Goettingen, Germany

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

resident in _____
City

to represent me/us at the Annual General Shareholders' Meeting of Sartorius AG convened for
April 7, 2016, 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 by midnight on **April 6, 2016 (24:00 hours Central European Time (CET) time of receipt)** to the following address: Sartorius Aktiengesellschaft c/o Better Orange IR & HV AG, Haidelweg 48, 81241 München, Germany, Fax: +49(0)89.889690633, 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 by midnight on **April 6, 2016 (24:00 hours CET time of receipt)** to the following address: Sartorius Aktiengesellschaft c/o Better Orange IR & HV AG, Haidelweg 48, 81241 München, Germany, Fax: +49(0)89.889690633, 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 AG