

Important Note:

The form provided below may be completed only by owners of ordinary shares as preference shares do not have any voting rights.

POWER OF ATTORNEY

Number of entrance card: _____

Number of ordinary shares: _____

First Name of Depositor

Last Name of Depositor

Depositor's City, Country and Postal Code/ZIP

Depositor's Street Address

IMPORTANT/URGENT MATERIAL: Please return by April 5, 2017

☐ Power of Attorney and Instructions to the Proxies Designated by the Company

I/we hereby grant power of attorney to each of the proxies designated by Sartorius Aktiengesellschaft, Ms. Marina Ciottariello and Ms. Leonie Helbling, both residents of Goettingen, Germany, and each entitled to grant substitute power of attorney, to represent me/us at the Annual General Shareholders' Meeting of Sartorius Aktiengesellschaft on April 6, 2017, by disclosing my/our name(s) and to exercise my/our voting right(s) according to my/our instructions given below.

or

In this case, please give instructions on how these proxies are to vote concerning the proposals of the management board listed in the items on the agenda below. If you do not place an X in any of the following boxes, the proxy(ies) will abstain from voting. For each item on the agenda, you may place an X **in only one of the boxes**. If both boxes are marked with an X, this will be considered invalid votes. (Please indicate your choice with an X)

☐ Voting by Mail

I/we will not be attending the Annual General Shareholders' Meeting of Sartorius Aktiengesellschaft in person and therefore vote by mail. For each item on the agenda, you may place an X **in only one of the boxes**. If both boxes are marked with an X, these will be considered invalid votes. (Please indicate your choice with an X)

(Please use a black or blue ink pen to complete and sign the form.)

Items on the Agenda

	YES	NO
2. Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft	<input type="checkbox"/>	<input type="checkbox"/>
3. Resolution on granting discharge to the members of the Executive Board for fiscal 2016	<input type="checkbox"/>	<input type="checkbox"/>
4. Resolution on granting discharge to the members of the Supervisory Board for fiscal 2016	<input type="checkbox"/>	<input type="checkbox"/>
5. Elections for the Supervisory Board		
5a. Election of Dr. Daniela Favoccia as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
5b. Election of Dr. Lothar Kappich as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
5c. Election of Ms. Ilke Hildegard Panzer as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
5d. Election of Prof. Dr. Dres. h.c. Arnold Picot as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
5e. Election of Prof. Dr. Thomas Scheper as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
5f. Election of Prof. Dr. Klaus Rüdiger Trützschler as a member of the Supervisory Board	<input type="checkbox"/>	<input type="checkbox"/>
6. Appointment of an auditor for fiscal 2017 as well as an auditor for the audit review of the first-half financial report of 2017	<input type="checkbox"/>	<input type="checkbox"/>

Place, Date

Signature of the person/s making a declaration acc. to §126b BGB (German Civil Code)
Please sign above

Please indicate your telephone number here in case we have any questions and need to contact you:

Phone number

Important Information

1. Power of attorney granted and voting instructions given to the proxies designated by the company

If you do not wish to attend the Annual General Shareholders' Meeting in person and have not granted a third party power of attorney, we offer you the opportunity to have the accountable proxies designated by the company to vote on your behalf.

For this purpose, please use the front of this form. If you would like to use this service, please send us the completed and signed form no later than on April 5, 2017, as the date of arrival:

- **by post** (regular mail) to:

Sartorius Aktiengesellschaft, c/o Better Orange IR & HV AG, Haidelweg 48, 81241 Munich, Germany

or

- **by fax** to: +49(0)89.889.690.633

or

- **by email** to: sartorius@better-orange.de

2. Voting by mail

Shareholders who do not wish to attend in person and do not wish to appoint a proxy may vote by mail. Please complete the voting-for-mail form on the front page. Voting by mail shall be in writing, in the form of text or in electronic form and must have been received by the company no later than the end of **April 5, 2017**. Please return this to the address, fax or email address stated in Section 1 "Power of attorney granted and voting instructions given to the proxies designated by the company" above.

Banks, institutes or companies that are equal to said banks by law (according to § 135, Subsection 10, and § 125, Subsection 5, of the German Stock Corporation Law [AktG]) or persons defined by § 135, Subsection 8, of AktG, in particular shareholders' unions, and have been granted power of attorney as proxies by the respective shareholder(s) may also vote by mail. Votes cast by mail on time may be changed or revoked up until the end of **April 5, 2017**, in writing, in the form of text or electronically using the address stated above. The date on which this mail arrives at the company's registered address shall determine whether said mail has been received on time.

Further information on attending the Annual General Shareholders' Meeting, appointment of proxies and on voting by mail

Forms for granting power of attorney or assigning a person to represent your rights (in particular voting rights, including granting power of attorney and issuing instructions to proxies designated by the company) are included on each entrance card. In addition, these forms are available on our Internet site at www.sartorius.com/hauptversammlung and can be ordered by mail, by fax or email at the address stated above.

If you grant power of attorney to more than one proxy, the company is entitled to reject one or several of these proxies.

Even after you have granted power of attorney to the proxies designated by the company, you are still entitled to attend the Annual General Shareholders' Meeting in person. If you as a shareholder attend the Annual General Shareholders' Meeting in person, this will end the appointment of the proxies designated by the company; in this case, the proxies appointed by the company will not exercise your respective shareholder's participation and voting rights. Registration in person by a third-party proxy at the registration desk in the Lokhalle, Goettingen, Germany, on the date of the Annual General Shareholders' Meeting on April 6, 2017, does not automatically revoke the power of attorney and instructions issued to the proxies designated by the company. In this case, you will need to revoke said power of attorney issued to these company-designated proxies in the form of text. For this purpose, please use the corresponding form, which you will find on the Internet site of the company at www.sartorius.com/hauptversammlung. As an alternative, you can obtain this form also at the registration desks in the Lokhalle in Goettingen, Germany, among other places.

If you have granted power of attorney to the proxies designated by Sartorius Aktiengesellschaft, but have not issued any instructions, these proxies cannot exercise your voting rights at the Annual General Shareholders' Meeting. Your votes will not be counted if you fail to issue any instructions or if your instructions are unclear.

If the proxies designated by the company are granted power of attorney, their names will be additionally recorded in the attendance list for the Annual General Shareholders' Meeting.

If you need to read the countermotions/election proposals of other shareholders required to be made available, you may do so by accessing our Internet site at www.sartorius.com/hauptversammlung. You can endorse a countermotion that is exclusively designated to reject a particular proposal presented by the management board, by giving instructions to vote against the management board's proposal.

Please note that when exercising voting rights, the proxies designated by Sartorius Aktiengesellschaft are exclusively bound by the instructions given to them. These proxies are not available for exercising other shareholder's rights, such as asking questions or presenting motions or addressing statements to the Annual General Shareholders' Meeting.

Furthermore, the proxies will not vote on wider-ranging countermotions or on any other motions submitted during the Annual General Shareholders' Meeting. In these cases, they will abstain from voting. If you wish to have your shareholder's rights exercised beyond the scope described above for proxies designated by the company, you will be required to vote in person or through a third-party proxy (not the proxies designated by the company) to whom you have granted power of attorney.