

Invitation

Sartorius Aktiengesellschaft
2020 Annual General Shareholders' Meeting
(Virtual General Meeting)

ISIN DE0007165607 and
ISIN DE0007165631

We cordially invite the shareholders of Sartorius Aktiengesellschaft to the Annual General Shareholders' Meeting on Friday, June 26, 2020, 10:00 a.m. (CEST), to be live-streamed from Otto-Brenner-Strasse 20, 37079, Göttingen, Germany, on the following terms:

Based on Article 2 of the German Act on Mitigating the Consequences of the Covid 19 Pandemic in Civil, Insolvency and Criminal Procedure Law dated March 27, 2020 (COVID 19 Act), the Annual General Shareholders' Meeting shall be held without the physical presence of shareholders or their proxies (with the exception of the voting proxies designated by the company) as a

virtual general meeting.

For the shareholders and their proxies (with the exception of the voting proxies designated by the company), there shall be no right or opportunity to attend the meeting on site.

I. Agenda

1. Presentation of the approved annual financial statements of Sartorius Aktiengesellschaft and the endorsed consolidated financial statements for the year ended December 31, 2019; the combined management report for Sartorius Aktiengesellschaft and the Group, together with the explanatory report, included therein, of the Executive Board concerning the disclosures according to § 289a, Subsection 1, and § 315a, Subsection 1, of the German Commercial Code (HGB); as well as together with the report of the Supervisory Board for fiscal 2019

The documents mentioned above are published on the Internet at the following address: www.sartorius.com/shareholders-meeting

2. Resolution on the appropriation of the retained profit of Sartorius Aktiengesellschaft

The Executive Board and the Supervisory Board will propose to appropriate the retained profit of €117,641,275.26 for fiscal 2019 as follows:

Payment of a dividend of €0.35	
per dividend-bearing ordinary share	= €11,974,278.40
Payment of a dividend of €0.36	
per dividend-bearing preference share	= €12,303,384.48
Balance of unappropriated profit carried forward to the new statement	
	€93,363,612.38
Total:	€117,641,275.26

Should the number of dividend-bearing shares change by the time the resolution on the appropriation of the retained profit is passed, a correspondingly adapted resolution proposal shall be submitted to a vote. Dividends will be paid out as of July 1, 2020.

3. Resolution on granting discharge to the members of the Executive Board for fiscal 2019

The Executive Board and the Supervisory Board will propose to grant discharge to the members of the Executive Board for fiscal 2019.

4. Resolution on granting discharge to the members of the Supervisory Board for fiscal 2019

The Executive Board and the Supervisory Board will propose to grant discharge to the members of the Supervisory Board for fiscal 2019.

5. Resolution on the amendment of the Articles of Association

The Executive Board and the Supervisory Board will propose to amend the Articles of Association as follows:

- Article 14, Subsection (1), of the Articles of Association will be amended as follows:
The Annual General Shareholders' Meeting shall be convoked by the Executive Board or the other persons legally authorized for this purpose.
- Article 14, Subsection (2), of the Articles of Association will be amended as follows:
The German legal provisions shall apply to the period of notice of convocation.
- Article 14, Subsection (3), of the Articles of Association will be deleted.

4. Article 15, Subsection (1), of the Articles of Association will be amended as follows:
Only those shareholders shall be entitled to attend the Annual General Shareholders' Meeting and to exercise their voting rights who have registered prior to the Annual General Shareholders' Meeting in accordance with the following provisions and have provided proof of their entitlement to participate.
5. Article 15, Subsection (2), of the Articles of Association shall be amended as follows:
Registration to attend the Annual General Shareholders' Meeting and proof of entitlement to participate shall be in text form in German or English (§ 126b of the German Civil Code BGB), unless a different form is permitted in the notice of convocation. For this purpose, proof of shareholding in accordance with § 67c, Subsection 3, of AktG shall be sufficient in any case in order to submit said proof of entitlement to participate. Proof of shareholding shall refer to the legally stipulated date before the Annual General Shareholders' Meeting (record date).
6. Article 15, Subsection (3), of the Articles of Association will be amended as follows:
Registration and proof shall be received by the company at the address specified for this purpose in the invitation within the respective legally stipulated period before the Annual General Shareholders' Meeting. The notice convening the meeting may also provide for shorter periods of notice, to be measured in days
7. Article 15, Subsection (4), of the Articles of Association will be amended as follows:
Voting rights may be exercised by a proxy. With regard to the form in which the power of attorney is granted, revoked and/or proven, the notice of convocation may provide for simplifications compared to the legally prescribed form; otherwise, the German statutory provisions shall apply. The provisions of § 135 of AktG shall remain unaffected thereby.
8. Article 15, Subsection (7), of the Articles of Association will be amended as follows:
Members of the Supervisory Board may participate in the Annual General Shareholders' Meeting by means of video and audio transmission if the respective member of the Supervisory Board is prevented for health, professional or personal reasons from attending the meeting in person at the designated meeting place.
9. The following clause shall be added as Article 19, Subsection (4), to the Articles of Association:
The Executive Board is authorized, in accordance with the provisions of § 59 of the German Stock Corporation Law (AktG), to pay an advance to the shareholders after the end of the fiscal year on the expected net retained profit for the year.

10. The Executive Board shall be instructed to file these amendments to the Articles of Association with the commercial register only on the cut-off date specified in § 26j, Subsection 4, of the Introductory Act to the German Stock Corporation Law (AktG) with regard to the applicability of § 67c of AktG (i.e., for registration on September 3, 2020).

6. Elections for the Supervisory Board

The term of Supervisory Board member Prof. David Ebsworth, PhD, shall expire at the end of this Annual General Shareholders' Meeting. Prof. David Ebsworth was appointed as a representative of the shareholders to the Supervisory Board effective January 1, 2020, for a term up to the end of the next Annual General Shareholders' Meeting by court resolution of the company's local court of registration Amtsgericht Göttingen to replace Dr. Guido Oelkers, who had resigned from his position as a member of the Supervisory Board with effect upon the expiration of December 31, 2019. For this reason, election of a Supervisory Board member is required to be held at the Annual General Shareholders' Meeting on June 26, 2020, pursuant to Article 8, Subsections (1) and (2), of the Articles of Association of Sartorius Aktiengesellschaft.

Pursuant to § 96, Subsection 1, and § 101, Subsection 1, of the German Stock Corporation Law (AktG) and to § 7, Subsection 1, sentence 1, item no. 1, of the German Codetermination Law (MitbestG), the Supervisory Board of Sartorius Aktiengesellschaft consists of six members representing the shareholders and of six members representing the employees, respectively, and, according to § 96, Subsection 2, sentence 1, of AktG, of at least 30% women (i.e., at least four) and of at least 30% men (i.e., at least four). The Supervisory Board must meet these minimum quotas on the whole, unless the Board's shareholder representatives or employee representatives object thereto pursuant to § 96, Subsection 2, sentence 3, of AktG. The shareholder representatives and the employee representatives each resolved to have the individual quotas separately met. Therefore, at least two women and at least two men must be appointed to the Supervisory Board for each of the sides constituted by the shareholder representatives and the employee representatives. These minimum quotas are already met regardless of the outcome of the election.

Said election shall be held in accordance with Article 8, Subsection (2), of the Articles of Association of Sartorius Aktiengesellschaft in a period not to exceed the end of the particular Annual General Shareholders' Meeting that passes a resolution on discharge of the Board members for the fourth fiscal year following the beginning of the term served on the Supervisory Board, with the fiscal year in which said term begins not counted.

The Annual General Shareholders' Meeting shall not be bound by the election proposals when electing members representing the shareholders to serve on the Supervisory Board.

At the recommendation of the Nomination Committee, the Supervisory Board proposes that Prof. David Ebsworth be elected as a representative of the shareholders to the Supervisory Board for a term up to the end of the Annual General Shareholders' Meeting that passes a resolution on granting discharge for fiscal 2021.

This election proposal is in line with the profile of skills and expertise for the Supervisory Board and with the objectives that this Board has set concerning its composition.

Prof. David Ebsworth, a resident of Overath, Germany, is currently engaged as a consultant for various healthcare companies and for financial investors.

He is a member of the following supervisory boards required to be set up under law or of comparable German national or international supervisory committees or oversight and control bodies of commercial enterprises:

- Verona Pharma plc, U.K. (*stock exchange listed*) – Chairman of the Board of Directors (*non-executive*)
- Pharma Investments SA, Luxembourg – Member of the board "Conseil d'Administration" (*non-executive*)
- Actimed Therapeutics Ltd, U.K. – Chairman of the Board of Directors (*non-executive*)

- Kyowa Kirin International plc, U.K. – member of the Board of Directors (*non-executive*)
- Opterion Health AG, Switzerland – Chairman of the Board "Verwaltungsrat" (*non-executive*)
- Interpharma Investments Ltd, British Virgin Islands – member of the Board of Directors (*non-executive*)

In the appraisal of the Supervisory Board, there are no material personal or business relationships within the meaning of the German Corporate Governance Code between Prof. David Ebsworth and Sartorius Aktiengesellschaft, its Group companies, the corporate bodies of Sartorius Aktiengesellschaft or a shareholder owning a material interest in Sartorius Aktiengesellschaft.

The curriculum vitae of Prof. David Ebsworth, which also lists his major activities in addition to his Supervisory Board mandate, is annexed to this invitation.

7. Appointment of an auditor for fiscal 2020 as well as an auditor for the audit review of the first-half financial report of 2020

Upon recommendation of the Audit Committee, the Supervisory Board will propose to appoint KPMG AG Wirtschaftsprüfungsgesellschaft, Hanover, Germany, as the auditors for fiscal 2020 and for the audit review of the first-half financial report of 2020.

II. Total Number of Shares and Voting Rights at the Time of Convocation

The company issued 74,880,000 bearer no-par shares, divided into 37,440,000 ordinary shares and 37,440,000 non-voting preference shares at the time the Annual General Shareholders' Meeting was convoked. At this time of convocation, the number of voting rights is 37,440,000.

The number entitled to participate is 68,388,292 shares as the company holds 3,227,776 ordinary shares and 3,263,932 preference shares; the company has no rights derived from these.

III. Further information on convocation

1. Virtual general meeting

Based on the COVID 19 Act, the Executive Board of the company with the approval of the Supervisory Board decided to hold the Annual General Shareholders' Meeting without the physical presence of the shareholders or their proxies (virtual general meeting).

The right of shareholders to follow this virtual general meeting on the Internet, as well as the rights of shareholders during this meeting, including the prerequisites for exercising such rights – either by the shareholders themselves or by a proxy they appoint – will be described in more detail under the following lettered items.

a) Registration for the virtual general meeting

Shareholders and their appointed proxies may exercise their rights stated in b) to e) only if they have duly registered for the virtual general meeting. The registration procedure is described in detail in Section III.2. of this invitation.

b) Video and audio transmission

For shareholders and their proxies, the entire virtual general meeting, including the responses given to questions as well as votes, will be live streamed by video and audio transmission on a password-protected Internet service platform at www.sartorius.com/shareholders-meeting. Shareholders and/or their proxies will be sent personal access data required to watch this live stream after they have duly registered for the virtual general meeting. Transmission of the virtual general meeting does not enable any participation in the Annual General Shareholders' Meeting within the meaning of § 118, Subsection 1, sentence 2, of AktG. Rather, the rights of shareholders and their proxies during the virtual general meeting are derived from the following subsections c) to e).

c) Exercise of voting rights

Exercise of voting rights by shareholders and their proxies shall be by way of voting by mail (postal voting) or by the voting proxies designated by the company who are duly instructed for this purpose by the shareholders. Both options for exercising voting rights are described in detail in Section III.3.

d) Opportunity to ask questions

Shareholders and their proxies can address their questions to the company administration about matters concerning the company insofar responses to these questions are required for appropriately assessing and completing an item on the agenda. To ensure that questions are answered under the difficult conditions of the COVID-19 pandemic, the Executive Board decided, with the approval of the

Supervisory Board, that for organizational reasons such questions shall be submitted to the company in German no later than by the end of **June 23, 2020 (midnight, 24:00 hours (CEST))** in accordance with Article 2, Section 1, Subsection (2), bullet number 2 of the COVID-19 Act via the password-protected Internet service platform at www.sartorius.com/shareholders-meeting. The date on which a question or questions arrive at the company shall determine whether said question(s) has(have) been received on time. Any questions submitted by other means or later shall not be considered.

Pursuant to Article 2, Section 1, Subsection (2), of the COVID-19 Act, the Executive Board is entitled to decide at its due and free discretion as to which questions it will answer ahead of the general meeting on the company's Internet site.

e) Submitting objections

Shareholders or their proxies who have exercised their voting rights according to Section III.1.c) are entitled to submit an objection to a resolution of the Annual General Shareholders' Meeting (virtual general meeting) for the notary public's minutes via the password-protected Internet service platform at www.sartorius.com/shareholders-meeting from the beginning to the end of the Annual General Shareholders' Meeting for the notary's minutes.

2. Prerequisites for exercising shareholders' rights; registration procedure; record date

Those shareholders – including only those ordinary shareholders who have registered by the end of **June 19, 2020 (24:00 hours [CEST])** at the latest at the address stated below are entitled to exercise the shareholder rights described in Subsections b) to e) of Section III.1. Their registration must be in text form (§ 126b of the German Civil Code [BGB]) and must be submitted in German or English.

Furthermore, the shareholders must submit proof of their entitlement to exercise their particular shareholder rights described in subsections b) to e) in Section III.1. For this purpose, shareholders shall be required to obtain proof of their shareholding, issued in the form of text, from the depositing or custodial bank, where said proof shall refer to the beginning of **June 5, 2020 (0:00 hours CEST, the so-called "record date")**, and must be received by the company no later than by the end of **June 19, 2020 (midnight, 24:00 hours CEST)**, at the address stated below. This proof shall be in the form of text (§ 126b of BGB) and be in German or English.

In relation to the company, only those persons shall be considered shareholders entitled to exercise their respective shareholder rights described in Subsections b) to e) in Section III.1. who have furnished proof of their shareholding. Entitlement shall be governed exclusively by their respective shareholding reported on the record date. This record date shall not impose a freeze on the sale of shareholdings. Even in the event that a shareholder sells his or her shareholding completely or partially after the record date, the shareholder's respective shares owned as of the record date shall exclusively govern his or her entitlement to exercise his or her shareholder rights described in Subsections b) to e) in Section III.1; i.e., any sales of shares after the record date shall not have any effect on a shareholder's entitlement to exercise his or her shareholder rights described in Subsections b) to e) in Section III.1. The same shall apply to all purchases of shares after the record date. Persons who do not yet own any shares as of the record date and do not become shareholders until afterwards shall not be entitled to exercise their particular shareholder rights described in Subsections b) to e) in Section III.1. The record date shall not have any significance concerning entitlement to receive dividends.

Registration and proof of shareholding shall be submitted to the following registration address:

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

or by fax: +49(0)89.889.690.633
or by email: sartorius@better-orange.de

Once a shareholder's registration and proof of his or her shareholding have been received will the respective shareholder be sent the access data for use of the password-protected Internet service platform. We ask shareholders to ensure that they register and send their proof of shareholding to the company well in advance.

3. Procedure for voting; voting by proxy

Owners of preference shares are not entitled to vote at the Annual General Shareholders' Meeting (virtual general meeting). The following explanations for representation by proxy(ies) therefore apply only to holders of ordinary shares.

a) Voting by mail (postal voting)

Shareholders duly registered can exercise their voting right by way of voting by mail. For this purpose, they can cast their vote up to **June 25, 2020, 6:00 p.m. (CEST)** using the following postal address, fax number or email 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

As an alternative, they may cast, change or revoke their vote **until voting begins on the date of the Annual General Shareholders Meeting on June 26, 2020**, via the password-protected Internet service platform at **www.sartorius.com/shareholders-meeting**. The date and time at which their vote is received shall govern whether they have met the respective deadline.

Those shareholders who wish to cast their vote by way of voting by mail are requested to use the password-protected Internet service platform at **www.sartorius.com/shareholders-meeting** or the form for voting by mail sent to them. As an alternative, the form for voting by mail will be sent to shareholders or their proxies at any time requested and can also be downloaded from the Internet at **www.sartorius.com/shareholders-meeting**

b) Voting by proxy

Shareholders who do not wish to personally exercise their voting rights may have these exercised by appointed proxies, such as a depositing or custodial bank, a union of shareholders, proxies designated by the company who shall be bound by the respective shareholders' specific instructions, or another person of their choice. In the event that a shareholder elects to be represented by a voting proxy, he or she shall still be required to register his or her shareholding by the deadline specified and submit proof thereof according to the requirements stated above. If a shareholder grants power of attorney to more than one proxy, the company is entitled to reject one or several of these proxies.

Granting of power of attorney to proxies or to the proxies designated by the company can be granted electronically using the password-protected Internet service platform at **www.sartorius.com/shareholders-meeting** or in the form of text (§ 126b of BGB).

The postal address, fax number and email address specified under Section III.3.a) and the password-protected Internet service platform at www.sartorius.com/shareholders-meeting will be available for granting power of attorney to the company and for transmitting proof of power of attorney granted to the proxy, as well as for voting by the proxy and for changing or revoking said power of attorney, in each case up to the deadlines specified there.

On the day of the virtual general meeting, powers of attorney may be electronically submitted, amended or revoked only by using the password-protected Internet service platform accessible at www.sartorius.com/shareholders-meeting up until the start of voting.

Proxies cannot attend the virtual general meeting either. They can exercise the voting rights for the shareholders they represent only within the scope of their respective power of attorney by way of a postal vote or by issuing substitute power of attorney to the company's designated proxies bound by the instructions of the respective shareholders.

Use of the password-protected Internet service platform by the proxy requires that the proxy receive the relevant access data, which will be sent to the shareholder after due registration for the virtual general meeting and proper proof of share ownership, from the person granting the power of attorney. If power of attorney is granted to banks, shareholders' unions, other intermediaries covered by § 135 of AktG or to persons and institutes equal to the same pursuant to § 135 of AktG, it is sufficient, however, if the declaration of power of attorney is held by the authorized proxy so that said power of attorney can be evidenced; in this case, the declaration of power of attorney must be complete and contain only the statements associated with the exercise of voting rights.

With regard to the revocation or amendment of a power of attorney, the foregoing statements with respect to granting of such power of attorney shall apply accordingly.

A form for power of attorney and further information on granting power of attorney, along with access data for using the password-protected Internet service platform, will be sent to duly registered persons. On request, the power of attorney form will be sent to shareholders or their proxies at any time and can furthermore be downloaded from www.sartorius.com/shareholders-meeting. Shareholders are requested to grant power of attorney by preferably submitting their completed power of attorney form using the password-protected Internet service platform at www.sartorius.com/shareholders-meeting or to use the form provided by the company for granting power of attorney.

Sartorius Aktiengesellschaft will offer holders of ordinary shares who have duly registered for the Annual General Shareholders' Meeting (virtual general meeting) the additional opportunity to appoint proxies designated by the company and bound by the instructions issued even before the Annual General Shareholders' Meeting takes place. In the event that proxies are appointed through power of attorney granted by the company, said proxies shall exercise the respective voting rights according to the binding instructions given by the respective shareholders. The proxies designated by the company will not be able to exercise a shareholder's voting rights if said shareholder has failed to issue any instructions. Power of attorney shall be granted and voting instructions shall be given in the form of text.

Those shareholders who wish to authorize and instruct the proxies designated by the company are requested to use the password-protected Internet service platform at www.sartorius.com/shareholders-meeting or the form sent to them for granting power of attorney and issuing instructions. Alternatively, the power of attorney and instruction form will also be sent to shareholders or their proxies at any time upon request and can also be downloaded from the company's Internet site at www.sartorius.com/shareholders-meeting.

Powers of attorney and instructions issued by duly registered shareholders to voting proxies designated by the company must be received by the company at the postal address, fax number or email address specified in Section III.3.a) or via the password-protected Internet service platform at www.sartorius.com/shareholders-meeting by the deadlines specified therein; the same shall apply to the amendment or revocation of power of attorney and of instructions. The date of receipt by the company shall determine whether said power of attorney and instructions issued have been received on time in each case.

Information on granting power of attorney and giving instructions to voting proxies designated by the company is also available to our shareholders on the company's Internet site at www.sartorius.com/shareholders-meeting.

The appointment of voting proxies to submit motions, ask questions and to file objections is excluded.

4. Further shareholders' rights

a) Right to put additional items on the agenda pursuant to § 122, Subsection 2, of AktG

Shareholders whose shares total one twentieth of the share capital or attain a proportionate amount of €500,000.00 may request that items be added to the agenda of the shareholders' meeting and be announced (§ 122, Subsection 2, sentence 1, in conjunction with Subsection 1, of AktG). Their request must be addressed in writing to the Executive Board of the company. Each such new item must be accompanied by a justification or a draft resolution.

Shareholders so requesting said addition shall submit proof according to § 122, Subsection 2, in conjunction with Subsection 1, of AktG that they have owned the company's shares for at least 90 days before the date of receipt of their request and that they will continue to hold said shares up until the Executive Board's decision on said request. For calculation of the number of days of this deadline, § 70 and § 121, Subsection 7, of AktG shall be observed.

Said request must be received by the company no later than by the end of **May 26, 2020 (midnight, 24:00 hours CEST)**, at the following address:

Sartorius Aktiengesellschaft
Executive Board
Otto-Brenner-Strasse 20
37079 Göttingen, Germany

b) Motions from shareholders and election proposals pursuant to § 126, Subsection 1, and § 127 of AktG

Motions from shareholders against a proposal of the Executive Board and the Supervisory Board concerning a particular item on the agenda in line with § 126, Subsection 1, of AktG, as well as against proposals of shareholders for the election of members to the Supervisory Board or for the appointment of auditors pursuant to § 127 of AktG must be directed to the following address only:

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

or by fax: +49(0)89.889.690.633
or by email: sartorius@better-orange.de

Motions sent to addresses other than the above will not be considered. Shareholders' motions to be made available concerning the agenda will be published, including the name(s) of the respective shareholder(s)

and a justification of this motion, without undue delay after its receipt, on the company's Internet site at www.sartorius.com/shareholders-meeting, provided that said motions together with a justification thereof are received by the company no later than by the end of **June 11, 2020 (midnight, 24:00 hours CEST)**. Opinions given by the company's management board concerning these motions will likewise be made available at this Internet address.

The company shall be entitled to refrain from publishing a countermotion along with its justification if one of the requirements set forth in § 126, Subsection 2, of AktG is met; requirements justifying the invocation of this right are met, for instance, if a countermotion will lead to the approval by a shareholders' meeting of a resolution that is against the law or contrary to the company's Articles of Association. Justification of a countermotion does not have to be published if the entire text exceeds 5,000 characters.

The statements made above apply to a shareholder's proposals for the election of members to the Supervisory Board or for the appointment of auditors, with the requirement that an election proposal does not have to be justified. Except for cases named in § 126, Subsection 2, of AktG, the company may also refrain from publishing proposals for election of members to the Supervisory Board and for appointment of the auditors if said proposals fail to list the name, job currently held and city of residence of the person(s) suggested, and, in the case of proposals for election of Supervisory Board members, if said proposals fail to include information on their membership in any legally constituted supervisory boards or advisory boards.

c) Right to receive information pursuant to § 131, Subsection 1, of AktG in conjunction with Article 2, Section 1, Subsection 2, of the COVID-19 Act

Shareholders and their proxies can address their questions to the company administration about matters concerning the company insofar as responses to these questions are required for appropriately assessing and completing an item on the agenda. To ensure questions are answered under the more difficult conditions of the COVID-19 pandemic, the Executive Board decided, with the approval of the Supervisory Board, that for organizational reasons such questions are to be submitted in German to the company, in accordance with Article 2, Section 1, Subsection 2, of the COVID-19 Act, no later than by the end of **June 23, 2020 (midnight, 24:00 hours (CEST))** in accordance with Article 2, Section 1, Subsection (2), bullet number 2 of the COVID-19 Act via the password-protected Internet service platform at www.sartorius.com/shareholders-meeting.

Pursuant to Article 2, Section 1, Subsection 2, of the COVID-19 Act, the Executive Board is entitled to decide at its due and free discretion as to which questions it answers and how. The Executive Board reserves the right to answer questions on the company's Internet site ahead of the virtual general meeting.

Explanations on the above-stated rights of shareholders in accordance with § 122, Subsection 2; § 126, Subsection 1; § 127; § 137; and § 131, Subsection 1; of AktG, in conjunction with Article 2, Section 1, Subsection 2, of the COVID-19 Act, where applicable, can also be downloaded from the company's Internet site at www.sartorius.com/shareholders-meeting.

5. Information on the company's Internet site (§ 124a of AktG)

Extensive information about the Annual General Shareholders' Meeting is given on the company's Internet site at www.sartorius.com/shareholders-meeting. This information includes, inter alia, the content of the convocation; an explanation if no resolution is to be adopted for a specific item on the agenda; the documents to be made available to the shareholders' meeting; the total number of shares and the voting rights at the time of convocation, including separate information on the total number of each class of share; the forms that are to be used for voting by proxy or by mail, unless these forms have already been sent directly to the shareholder(s); and proof of publication of the invitation in the German Federal Gazette (Bundesanzeiger).

6. Information on data protection

a) Legal basis for processing personal data

Within the scope of conducting the Annual General Shareholders' Meeting, the company processes the personal data of its shareholders and those of their proxy(ies), if any (in particular the name, address, email address, number of shares, class of share, type of share ownership and the ID number of confirmation of registration) on the basis of the data protection laws in force in order to enable shareholders to exercise their rights pursuant to AktG and the Articles of Association of Sartorius Aktiengesellschaft within the context of the Annual General Shareholders' Meeting. Processing of personal data is mandatory in order to properly prepare for, conduct and follow up on the Annual General Shareholders' Meeting and to enable shareholders to participate in the Annual General Shareholders' Meeting in accordance with Sections 118 et seq. of AktG. The responsible entity for processing said personal data is Sartorius Aktiengesellschaft (Otto-Brenner-Strasse 20, 37079 Göttingen, Germany) within the meaning of Art. 4, no. 7, of the European General Data Protection Regulation ("GDPR"). The legal basis for processing

personal data is Article 6(1), sentence 1, letter (c), of the European General Data Protection Regulation (GDPR) in conjunction with § 67 and §118 et seq. of AktG, as well as in conjunction with Section 1 of the COVID-19 Act. In addition, data processing that is useful for the organization of the virtual general meeting may be carried out on the basis of overriding legitimate interests (Art. 6 (1), sentence 1, letter (f) of GDPR). As a rule, the company receives the shareholders' personal data via the registration office from the intermediary that the shareholders have appointed to hold their shares in custody (usually the custodial bank).

b) Transfer of personal data

The company engages various service providers for the purpose of holding the Annual General Shareholders' Meeting. The company provides them with said personal data only on a need-to-know basis in order for them to provide the requested service. The service providers are obligated to process said data solely in accordance with the company's instructions. In addition, personal data is made available to the shareholders and shareholder proxies in connection with the Annual General Shareholders' Meeting, by name in the list of participants, within the framework of the statutory regulations.

Moreover, the company transfers personal data of shareholders and of their proxies, if any, to further recipients outside the company, with said recipients processing these data under their own responsibility, Art. 4, no. 7 of GDPR. Such recipients may be, for example, public authorities due to legal requirements.

c) Storage of personal data

Personal data are stored as long as this is required by law or Sartorius Aktiengesellschaft has a legitimate interest in said storage, for example, any interest resulting from liability risks arising from applicable law. Afterwards, personal data are erased.

The period of storage for personal data collected in connection with the Annual General Shareholders' Meeting and in compliance with the legal obligations regarding provision of evidence and retention (e.g., in AktG, HGB, German Tax Code) is regularly up to three years. Data stored in the share register must be regularly stored for ten years after any sale of shares.

d) Rights regarding personal data

Data subjects shall have the following rights at any and all times concerning the processing of their personal data: right of information and access of personal data, rectification of said personal data, right to restriction of processing, right to object and the right to erasure ("right to be forgotten") in accordance with Chapter 3 of GDPR. These rights can be claimed from Sartorius Aktiengesellschaft free of charge through said company's

Data Protection Officer using the contact details in Section III.7 e) of this Invitation to the Annual General Shareholders' Meeting.

Furthermore, data subjects shall have the right to lodge a complaint with a data protection supervisory authority in accordance with Art. 77 of GDPR.

e) Data Protection Officer

The Data Protection Officer of the company shall be available to data subjects according to the following contact details in order to serve as a contact for matters related to data protection:

Sartorius Corporate Administration GmbH
Datenschutzbeauftragter (Data Protection Officer)
Otto-Brenner-Strasse 20
37079 Göttingen, Germany

or by email: datenschutz@sartorius.com

Göttingen, May 2020

Sartorius Aktiengesellschaft
The Executive Board

Annex to I.6 Elections for the Supervisory Board

Curriculum Vitae

Prof. David Ebsworth, PhD, resident of Overath, Germany
Born in 1954

Professional Profile:

Since 2015: Provision of consultancy services to selected healthcare companies and financial investors, as well as angel investor in selected healthcare companies

Since 2018:
Member of the Board "Conseil d'administration" (*non-executive*) of Pharma Investments SA, Luxembourg

Since 2018:
Chairman of the Board of Directors (*non-executive*) of Actimed Therapeutics Ltd, U.K.

Since 2017:
Member of the Board of Directors (*non-executive*) of Kyowa Kirin International plc, U.K.

Since 2016:
Chairman of the Board "Verwaltungsrat" (*non-executive*) of Opterion Health AG, Switzerland

Since 2016:
Member of the Board of Directors (*non-executive*) of Interpharma Investments Ltd, British Virgin Islands

Since 2016:
Visiting Professor, Business School of the University of Surrey, Guildford, U.K.

Since 2014: Chairman of the Board of Directors (*non-executive*) of Verona Pharma plc, U.K. (*stock market listed*)

2009–2015: Galenica AG, Switzerland

2014–2015:
Project support for CEO of Vifor Pharma and CEO of Galenica Santé until retirement

2012–2014:
CEO and Chairman of the Corporate Executive Committee (Zurich and Bern, Switzerland)

2009–2014:
CEO of Vifor Pharma AG and member of the Corporate Executive Committee of Galenica AG

2003–2009: Various positions as non-executive director and consultant to the industry and financial investors

2002–2003: CEO of Oxford Glycosciences Ltd, U.K.

1983–2002: Bayer AG, Germany

2000–2002:
President and General Manager, Pharmaceuticals

1995–1999:
President, North American Pharmaceutical Division

1983–1995:
Various positions (focal areas: Marketing, Sales, Business Operations)

Educational Background:

- B.Sc. in Chemistry and German (1976), University of Surrey, Guildford, U.K.
- Ph.D. in Comparative Industrial Relations (1980), University of Surrey, Guildford, U.K.
- Doctor of Humane Letters (2000), Honorary Degree, University of New Haven, Connecticut, USA

Major Activities Beyond Supervisory Board Position:

- Consultant for various healthcare companies and financial investors
- Visiting Professor, Business School of the University of Surrey, Guildford, U.K.
- Positions held within the meaning of § 125, Subsection 1, sentence 5, of the German Stock Corporation Law (AktG)

Positions Held Within the Meaning of § 125, Subsection 1, Sentence 5, of the German Stock Corporation Law (AktG):

- Verona Pharma plc, Großbritannien (*börsennotiert*) – Chairman of the Board of Directors (*non-executive*)
- Pharma Investments SA, Luxembourg – member of the board “Conseil d’administration” (*non-executive*)
- Actimed Therapeutics Ltd, U.K. – Chairman of the Board of Directors (*non-executive*)
- Kyowa Kirin International plc, U.K. – member of the Board of Directors (*non-executive*)
- Opterion Health AG, Switzerland – Chairman of the Board “Verwaltungsrat” (*non-executive*)
- Interpharma Investments Ltd, British Virgin Islands – member of the Board of Directors (*non-executive*)

■ Sartorius Annual Report Available as Both a Download Version and an Online Version

Sartorius is increasingly using online formats for its communication and thus offers you, as a user, additional functions, such as a full-text search and a download center for tables.

Moreover, the Sartorius Annual Report is available for you to download as a PDF file from our website at www.sartorius.com/shareholders-meeting.

- Sartorius relaunched its brand as of February 17, 2020, and in this context also introduced a new visual look. For this reason, the present invitation to the 2020 Annual General Shareholders' Meeting is issued in the new layout, while our annual report is published in our former layout as the editorial deadline was before the relaunch date on the February 17.

Sartorius AG
Otto-Brenner-Straße 20
37079 Göttingen

Telefon 0551 308 0
Fax 0551 308 3955

www.sartorius.com