

Total Number of Shares and Voting Rights at the Time of Convocation – Information According to § 124a, No. 5, of AktG

On the date that the announcement of the convening of the Annual General Meeting was published in the Federal Gazette (Bundesanzeiger), the company had issued 74,880,000 no-par value bearer shares, divided into 37,440,000 ordinary shares and 37,440,000 non-voting preference shares. Preference shares do not confer voting rights except in cases specified by law. To the extent that shareholders have voting rights, each ordinary or preference share entitles the holder to one vote at the Annual General Meeting.

Unless the agenda of this Annual General Meeting is amended via a separate announcement to include additional items on which preference shares are entitled to vote, only ordinary shares are entitled to vote at this Annual General Meeting. The total number of voting rights therefore corresponds to the total number of ordinary shares and amounts to 37,440,000 on the date that the announcement of the convening of the Annual General Meeting was published in the Federal Gazette (Bundesanzeiger).

Pursuant to Section 71b of the German Stock Corporation Act (AktG), the company does not have any voting rights from directly or indirectly held treasury shares. On the date that the announcement of the convening of the Annual General Meeting was published in Germany's Federal Gazette (Bundesanzeiger), the company held 3,201,331 ordinary shares and 2,626,379 preference shares in the company as treasury shares.

Sartorius Aktiengesellschaft