

# Total number of Shares and Voting Rights at the Time of Convocation – Information According to § 124a, No. 4, of AktG

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 68,415,862 shares as the company holds 3,213,991 ordinary shares and 3,250,147 preference shares; the company has no rights derived from these.

Sartorius Aktiengesellschaft