

TOTAL NUMBER of Shares and Voting Rights at the Time of Convocation

The company issued 18,720,000 bearer no-par shares, divided into 9,360,000 ordinary shares and 9,360,000 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 9,360,000. The number entitled to participate is 17,047,073 shares as the company holds 831,944 ordinary shares and 840,983 preference shares; the company has no rights derived from these.

Sartorius AG