

Explanation According to § 124a, No. 2, of AktG

At the meeting on February 13, 2020, the Supervisory Board endorsed the annual financial statements prepared by the Executive Board for the year ended December 31, 2019, and the consolidated financial statements for the year ended December 31, 2019; the annual financial statements are thus approved pursuant to § 172 of the German Stock Corporation Law (AktG). For this reason, it is not necessary pursuant to § 173 of AktG for the Annual General Shareholders' Meeting to approve the annual financial statements or to endorse the consolidated financial statements so that no resolution will be adopted for item no. 1 on the agenda.

Sartorius Aktiengesellschaft