

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

The Supervisory Board approved the annual financial statements and the consolidated annual financial statements prepared by the Executive Board; the annual financial statements are thus adopted. In this case, the law does not provide for the adoption of the annual financial statements or the approval of the consolidated financial statements by the Annual General Meeting. Instead, the aforementioned documents are to be made accessible to the Annual General Meeting in accordance with the legal regulations (Section 176(1)(1) of the German Stock Corporation Act, "Aktiengesetz," abbreviated as "AktG" in German). Accordingly, no resolution will be passed by the Annual General Meeting on agenda item 1.

Sartorius Aktiengesellschaft