

## Declaration of the Executive Board and of the Supervisory Board of Sartorius AG Concerning the Recommendations of the Government Commission on the German Corporate Governance Codex Pursuant to Section 161 of the German Stock Corporation Law ("Aktiengesetz")

1. The Executive Board and the Supervisory Board declare that Sartorius AG complied, and will continue to comply in the future, with the recommendations made by the Government Commission on the German Corporate Governance Code (GCGC) in the version as amended on December 16, 2019, and published by the German Federal Ministry of Justice and Consumer Protection in the official section of the German Federal Gazette ("Bundesanzeiger") since their publication, with the following exception:

In divergence from the newly introduced recommendations pursuant to G.10 of GCGC, the variable remuneration of the members of the Executive Board consists only to a minor extent of share-based remuneration components; furthermore, portions of the long-term remuneration with a multi-year assessment basis are available to the members of the Executive Board before the end of four years. The Supervisory Board believes that the terms of variable remuneration in place so far also provide for an incentive structure that is geared towards the sustainable and long-term development of the company. The Supervisory Board will observe and evaluate the implementation of the recommendations according to G.10 of GCGC on the market and, if necessary, propose changes to the remuneration system to the Annual General Shareholders' Meeting at a later time.

2. The Executive Board and the Supervisory Board further declare that Sartorius AG fully complied with the recommendations made by the Government Commission on the German Corporate Governance Code in the version as amended on February 7, 2017, and published by the German Federal Ministry of Justice and Consumer Protection in the official section of the German Federal Gazette ("Bundesanzeiger") on April 24, 2017, and May 19, 2017, respectively, throughout the period from the issuance of last year's Declaration of Compliance in December 2019 to the publication of the new version of the GCGC on March 20, 2020.

Göttingen, December 3, 2020

For the Supervisory Board



Dr. Lothar Kappich

For the Executive Board



Dr. Joachim Kreuzburg