

**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 Code Pursuant to Section 161 of the German Stock
Corporation Act ("Aktiengesetz"):**

The Executive Board and the Supervisory Board declare that Sartorius AG complied in the period since last year's Declaration of Compliance was issued on December 6, 2024, and will continue to comply in the future, with the recommendations made by the Government Commission on the German Corporate Governance Code (GCGC) as of April 28, 2022, as published by the Federal Ministry of Justice in the official section of the Federal Gazette on June 27, 2022, with the following exception:

In divergence from the Recommendation pursuant to G.10, sentence 1, of the GCGC, the variable remuneration of the members of the Executive Board – up to and including June 30, 2025, with the exception of the variable remuneration of the then Chairman of the Executive Board – consists only to a non-predominant extent of share-based remuneration components. The Supervisory Board believes that the existing structure of the variable remuneration, which corresponds to the remuneration system approved by the Annual General Meeting, also achieves an incentive structure geared to the sustainable and long-term development of the Company.

Goettingen, December 4, 2025

For the Supervisory Board



Dr. Lothar Kappich

For the Executive Board



Dr. Michael Grosse