

**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 § 161 of the German Stock Corporation Law ("Aktiengesetz")**

The Executive Board and the Supervisory Board declare that full compliance will be achieved with the recommendations made by the Government Commission on the German Corporate Governance Code and published by the German Federal Ministry of Justice in the official section of the German Federal Gazette ("Bundesanzeiger"), as amended on February 7, 2017.

Since last year's Declaration of Compliance was issued, Sartorius AG has complied with the recommendations in the valid version of the Government Commission on the German Corporate Governance Code to the full extent.

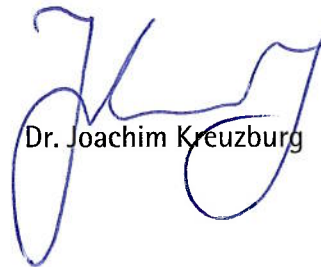
Goettingen, December 7, 2017

For the Supervisory Board



Dr. Lothar Kappich

For the Executive Board



Dr. Joachim Kreuzburg