

**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 electronic German Federal Gazette ("Elektronischer Bundesanzeiger"), as amended on May 13, 2013.

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.

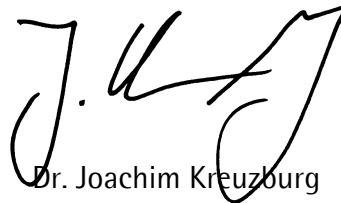
Goettingen, December 12, 2013

For the Supervisory Board



Prof. Dr. Dres. h.c. Arnold Picot

For the Executive Board



Dr. Joachim Kreuzburg

You will find the complete and up-to-date German Corporate Governance Code on the following web page in both English and German: www.corporate-governance-code.de; for this version in German only, visit www.ebundesanzeiger.de.