

**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 compliance 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 June 6, 2008.

However, this compliance applies with the restriction that a D&O insurance policy without any deductible by officers exists for the members of the Executive Board and of the Supervisory Board (Code Item 3.8, Paragraph 2) and that the remuneration of the Supervisory Board members is not individualized, but rather reported as a total subdivided into fixed and variable components, and as payments for services personally rendered, in the Notes to the Financial Statements and in the Corporate Governance Report (Code Item 5.4.6, Paragraph 3).

Since the issue of last year's Declaration of Compliance, the recommendations of the Government Commission of the German Corporate Governance Code have been complied with as specifically amended, subject to the restrictions already stated above.

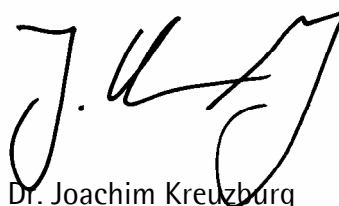
Goettingen, December 11, 2008

For the Supervisory Board



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

For the Executive Board



Dr. Joachim Kreuzburg

You can find a complete and up-to-date version of the German Corporate Governance Code on the following pages on the Internet: [www.ebundesanzeiger.de](http://www.ebundesanzeiger.de) and [www.corporate-governance-code.de](http://www.corporate-governance-code.de).