

Ad hoc announcement

according to Section 15, Subsection 1,
of the German Securities Trading Act (WpHG)



Sartorius plans to raise share capital by use of retained earnings / 4:1 "stock split"

Goettingen, February 25, 2016 - The Supervisory Board and the Executive Board of Sartorius AG today resolved to submit a proposal to the Annual General Shareholders' Meeting scheduled on April 7, 2016, to increase share capital by use of retained earnings for carrying out a so-called stock split for both ordinary and preference shares at a ratio of 4:1. For this purpose, share capital would be increased accordingly by the conversion of capital reserves. Each ordinary shareholder would receive three additional ordinary shares for every Sartorius ordinary share held, without an additional payment. Every preference share owner would likewise receive three additional preference shares for every Sartorius preference share held, without an additional payment.

This move will not have any impact on the amount of the company's equity as it will entail only a reclassification within equity capital. If the proposal is approved by a corresponding resolution passed by the Annual Shareholders' Meeting, the share capital of Sartorius AG would quadruple from 18,720,000 euros to 74,880,000 euros.

As part of this move to increase share capital, the share in dividends (surplus and minimum dividends) would also be adjusted for preference share owners. The holders of non-voting preference shares would receive an increased dividend (surplus dividend) of 0.01 euro per preference share from the distributable annual profit; however, the dividend would at least amount to 0.02 euro per preference share (minimum dividend).

This ad hoc announcement contains statements about the future development of the Sartorius Group. The content of these statements cannot be guaranteed as they are based on assumptions and estimates that harbor certain risks and uncertainties. This is a translation of the original German-language ad hoc announcement. Sartorius shall not assume any liability for the correctness of this translation. The original German ad hoc announcement is the legally binding version. Furthermore, Sartorius reserves the right not to be responsible for the topicality, correctness, completeness or quality of the information provided. Liability claims regarding damage caused by the use of any information provided, including any kind of information which is incomplete or incorrect, will therefore be rejected.

Contact

Petra Kirchhoff, Vice President of Corporate Communications
+49 (0)551.308.1686 | petra.kirchhoff@sartorius.com

Andreas Theisen, Director Investor Relations
+49 (0)551.308.1668 | andreas.theisen@sartorius.com