

Disclosure of inside information according to Article 17 MAR

Michael Grosse to become Chief Executive Officer of Sartorius AG

At its meeting today, the Supervisory Board of Sartorius AG resolved to appoint Dr. Michael Grosse as member of the Executive Board and Chief Executive Officer of the company with effect from July 1, 2025, for a term of three years. Dr. Joachim Kreuzburg, CEO since more than 20 years, had announced in July 2024 that he will not seek a further extension of his contract, which runs until November 2025, and will continue to lead Sartorius until mid-2025.

This notification according to Article 17 MAR 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 announcement. The Sartorius Group shall not assume any liability for the correctness of this translation. The original German ad hoc announcement is the legally binding version. Furthermore, the Sartorius Group 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.

Göttingen, December 6, 2024

Contact

Petra Kirchhoff

Head of Corporate Communications and Investor Relations

+49 (0)551.308.1686

petra.kirchhoff@sartorius.com

sartorius.com