

Sartorius appoints Gerry MacKay as Executive Vice President of Marketing, Sales and Services for its Laboratory Division

- **Experienced Biopharma manager to join Group Executive Committee**

Goettingen, Germany, March 9, 2017 – Sartorius, a leading international pharmaceutical and laboratory equipment provider, today announces the appointment of Gerry MacKay as Executive Vice President of Marketing, Sales and Services for its Lab Products & Services Division (LPS) with immediate effect. MacKay will report to CEO Dr. Joachim Kreuzburg, and be a member of the Group Executive Committee, the key management body of the Sartorius Group.

Gerry MacKay, who brings in a broad international background in the biopharma business, joined Sartorius in 2015 through the acquisition of BioOutsource, a contract testing services company, where he had served as CEO. Prior to his tenure with BioOutsource, he held senior level positions with Entegris, Mykrolis and Millipore.

"We are very pleased that Gerry has accepted this new challenging position at Sartorius. His considerable experience in the global biopharma market will be instrumental to strengthen the momentum on our LPS division's new strategic direction towards Biopharma and respective applications and technologies", said Sartorius CEO Joachim Kreuzburg.

MacKay assumes the role of his predecessor, Michael Melingo.



Image files

Gerry MacKay, Member of the Sartorius Group Executive Committee:

A Profile of Sartorius

The Sartorius Group is a leading international pharmaceutical and laboratory equipment provider with two divisions: Bioprocess Solutions and Lab Products & Services. Bioprocess Solutions with its broad product portfolio focusing on single-use solutions helps customers produce biotech medications and vaccines safely and efficiently. Lab Products & Services, with its premium laboratory instruments, consumables and services, concentrates on serving the needs of laboratories performing research and quality assurance at pharma and biopharma companies and on those of academic research institutes. Founded in 1870, the company earned sales revenue of more than 1.3 billion euros in 2016. 7,000 people work at the Group's 50 manufacturing and sales sites, serving customers around the globe.

Contact

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