

Chairman of the Supervisory Board Dr. Lothar Kappich Appointed Executor of Horst Sartorius' Will

Goettingen, September 25, 2017 – The local German court Goettingen (Amtsgericht) confirmed the appointment of Dr. Lothar Kappich as executor of Horst Sartorius' will. This new appointment became necessary after the former executor, Professor Arnold Picot, Ph.D., had passed away on July 9, 2017. This executorship with a period of administration defined up to 2028 remains unchanged in undivided form and covers around 50.1 percent of the ordinary shares of Sartorius AG.

Lothar Kappich, who holds a university degree in economics, has been a member of the Sartorius Supervisory Board since 2007 and, likewise as successor to Prof. Picot, Chairman of this board since July 19, 2017.

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. More than 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