

Sartorius Opens New Subsidiary in Peru

Lima, March 17, 2016 – Today, Sartorius, a leading international laboratory and pharmaceutical equipment supplier, officially opened its new sales office in Lima, Peru. Among the around 100 guests were representatives from local government, customers and business partners. The new subsidiary focuses on the expanding pharmaceutical market. It will offer a broad portfolio of single-use products and equipment used in bioprocessing. The office starts with four employees, additional hires are planned. In the past, Sartorius attended the Peruvian market through its sales site in Argentina.

In his opening remarks, Robert Eberlein, Managing Director of Sartorius Group Latin America, acknowledged the strategic importance of the new office: "We are very confident that Peru is the right place to invest in. From this new home base we will not only be able to serve the needs of the growing Peruvian pharma market better but also to expand our activities on the South American Pacific coast."

Photo



The Sartorius team in Lima

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 1.1 billion euros in 2015. Around 6,200 people work for the Group, which has its own manufacturing and sales sites in around 110 countries.

Contact

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