

Göttingen, Germany | September 18, 2025

Sartorius expands production of key components for cell and gene therapies in France

- New GMP facility for transfection reagents of French sub-group Sartorius Stedim Biotech in Illkirch
- Facility equipped to support next-generation technologies in cell and gene therapies

The life science group Sartorius has expanded the site of its French sub-group Sartorius Stedim Biotech in Illkirch, near Strasbourg. The state-of-the-art facility adds pharmaceutical-grade (GMP) production of transfection reagents, which are critical components for viral vector manufacturing for novel therapies. The investment underscores Sartorius' dedication to supporting customers in developing and producing cell and gene therapies, a rapidly growing segment in biopharmaceuticals.

"The expanded Illkirch site is a key manufacturing hub in our global network and enables us to even better help customers deliver advanced therapies to patients faster and more efficiently," said René Fáber, Member of the Sartorius Executive Board and CEO of Sartorius Stedim Biotech.

In total, the site has more than doubled in size and now covers more than 8,000 square meters of production, lab, and office space. The fully digitalized plant is designed to operate under the EU guidelines for sterile drug manufacturing (GMP Annex 1 standard) and combines flexibility for future growth with high environmental standards.

The Illkirch site expansion follows the recent completion of a major capacity upgrade in Aubagne, France. Sartorius is investing strategically in the expansion of its global development and manufacturing network in order to prepare the company for further organic growth and to ensure high delivery reliability and consistent product quality across all regions.

For more information, read the [full media release of Sartorius Stedim Biotech](#)
Click here for [images of Sartorius Stedim Biotech Illkirch](#)

A profile of Sartorius

Sartorius is a leading international partner to life sciences research and the biopharmaceutical industry. With innovative laboratory instruments and consumables, the Group's Lab Products & Services Division focuses on laboratories performing research and quality control at pharmaceutical and biopharmaceutical companies as well as academic research institutes. The Bioprocess Solutions Division, with its broad product portfolio focusing on single-use solutions, helps customers manufacture biotech medications, vaccines, and cell and gene therapies more safely, rapidly, and sustainably. Based in Göttingen, Germany, the company has a strong global reach with around 60 production and sales sites worldwide. Sartorius

regularly expands its portfolio through the acquisition of complementary technologies. In 2024, the company generated sales revenue of around 3.4 billion euros. More than 13,500 employees work for customers around the globe.

Visit our [Newsroom](#) or follow us on [LinkedIn](#).

Contact

Verena Sattel

External Communications Manager

+49 551 308 9261

verena.sattel@sartorius.com