

Göttingen, Germany | March 16, 2026

## Sartorius launches next-generation platform to boost efficiency in cell therapy production

- Eveo Cell Therapy Platform enables automated, multi-parallel intensified production of autologous cell therapies
- Compact design can increase output up to fourfold in existing cleanroom space
- 90 percent reduction in manufacturing costs possible in CAR-T processes
- Successful pilot project with CDMO ElevateBio proves applicability

The life science group Sartorius is launching an integrated system for the production and quality control (QC) of autologous cell therapies. The modular Eveo Cell Therapy Platform is designed to address structural manufacturing bottlenecks that limit scalability and patient access to transformative treatments such as CAR-T therapies. With its latest innovation, Sartorius introduces a fundamentally new approach to cell therapy production that strengthens the company's position in the dynamic advanced therapies market.

"Cell therapies are redefining medical possibilities. Yet, they remain accessible to only a small share of patients because manufacturing is complex, capacity-constrained and therefore expensive. With our integrated platform, we are tackling these structural bottlenecks head-on," said René Fáber, member of the Executive Board at Sartorius. "By combining raw materials, production equipment, software and QC assays into a scalable system, we enable our customers to expand output, significantly reduce cost per dose and shorten vein-to-vein time for patients."

Conventional manufacturing models for autologous cell therapies rely on centralized, purpose-built facilities, that require significant greenfield investment, staff recruitment, training and multi-year build-up phases. Nevertheless, around 60 percent of regulatory disruptions in cell therapy production are caused by manufacturing and process-related issues.

The new Eveo platform addresses these challenges by integrating cell selection, activation, gene modification, expansion, wash and concentration and final formulation into a closed system, using a variety of Sartorius solutions from critical raw materials to bioreactors and separation technologies. The system's compact design supports implementation in existing facilities with existing staff, allows for operation in lower-classification environments and supports centralized and decentralized manufacturing models.

Sartorius has collaborated with ElevateBio, a leading cell therapy CDMO, amongst other partners, to test out the Eveo platform. In a successful pilot, the system performed in line with design expectations, demonstrating its real-world applicability. Due to its multi-parallel configuration, a single operator can process eight patient batches at once in the same cleanroom space traditionally needed for two batches.

Assuming a manufacturing time of around seven days per batch and 50 production weeks per year, Eveo enables production of more than 350 doses per year in a footprint that today yields roughly 100 doses – an almost fourfold increase. Financial modelling confirms that the platform can reduce manufacturing costs by around 90 percent.

Quality control also represents a major bottleneck in cell therapy production and release. With Sartorius' capabilities in live cell imaging and high throughput flow cytometry included in the Eveo platform, the technology addresses key analytical challenges and enables accessibility.

Sartorius plans to begin taking orders for the Eveo Cell Therapy Platform in September 2026, with first deliveries expected in 2027. ElevateBio is a preferred partner and will be granted access to one of the first Eveo platforms for use with its manufacturing customers.

### **More information about Eveo Cell Therapy Platform**

[www.sartorius.com/en/pr/eveo-cell-therapy-platform](http://www.sartorius.com/en/pr/eveo-cell-therapy-platform)

### **A profile of Sartorius**

Sartorius is a leading international partner to the biopharmaceutical research and manufacturing industries. The Lab Products & Services Division focuses on innovative laboratory instruments and consumables for research and quality assurance laboratories in pharmaceutical and biopharmaceutical companies as well as academic research institutions. The Bioprocess Solutions Division supports customers with a broad product portfolio focused on single-use solutions for the safer, faster, and more sustainable production of biotech drugs, vaccines, and cell and gene therapies. With around 60 production and sales locations worldwide, the Göttingen-based company has a strong global presence. Sartorius regularly supplements its portfolio with acquisitions of complementary technologies. In 2025, the company generated sales revenue of around 3.5 billion euros. More than 14,000 employees serve customers around the globe.

Visit our [newsroom](#) and follow us on [LinkedIn](#).

### **Contact**

Leona Malorny  
Head of External Communications  
+49 551 308 4067  
[leona.malorny@sartorius.com](mailto:leona.malorny@sartorius.com)