

Göttingen, Germany | September 11, 2024

## Sartorius and G-CON collaborate to deliver sustainable, flexible cGMP facilities faster

- The collaboration presents G-CON POD® cleanrooms outfitted with fully configured and manufacturing-ready Sartorius ProcessGo® Solutions
- The collaboration increases speed and flexibility of new facility builds or site expansions to deliver life-saving therapies more sustainably and faster

The life science group Sartorius and modular cleanroom manufacturer G-CON are launching a collaboration to enable rapid deployment of manufacturing solutions for therapeutic drug production around the globe.

Sartorius will provide ready-to-order, configurable and modality-specific Sartorius ProcessGo® Kits inside G-CON PODs. G-CON is the industry leader in prefabricated modular cleanroom solutions, and its POD systems play a major role in drug production due to their readily deployable, flexible and scalable approach to cleanrooms.

The manufacturing-ready Sartorius ProcessGo® solutions include all the equipment and consumables necessary to execute a GMP-compliant production batch. These solutions are supported by Sartorius' automation platform, data acquisition and analytic software to realize peak operating efficiency. Sartorius ProcessGo® solutions are engineered to optimize space and maximize yield. When integrated into G-CON cleanroom PODs, sustainable manufacturing practices are promoted by reducing the facility's carbon footprint, largely through optimization of HVAC systems and efficiency of facility construction and commissioning. The consumables of the ProcessGo® Solution come from a library of standard parts, ensuring supply chain security and enhancing carbon efficiency by simplifying logistics.

"This collaboration with G-CON, a leader in modular cleanroom solutions, represents a significant step forward in our commitment to helping our clients meet their challenges in terms of speed, efficiency, quality and sustainability," said Jake Slivka, Head of Integrated Solutions Market Development at Sartorius. "By combining Sartorius' expertise in single-use bioprocessing with G-CON's state-of-the-art cleanroom technology, we aim to accelerate the development and manufacturing of life-saving therapies, ultimately leading to better health for more people."

"We are very excited to be working with Sartorius, a true industry leader in bioprocessing. The value of these collaborations continues to increase as our industry advances and adopts turnkey platform solutions for new manufacturing infrastructure," said Dennis Powers, SVP of Product & Strategy at G-CON. "The benefits of integrating innovative bioprocess technology platforms such as Sartorius ProcessGo® within G-

CON's modular turnkey cleanroom solutions have been well recognized and we look forward to advancing that even further in our work with Sartorius."

In time, this innovation will advance the industry by enabling and increasing accessibility to healthcare, environmental sustainability, global biopharmaceutical capacity, and localized manufacturing in developing regions.

**Discover the latest insights and technology in process intensification and attend a joint presentation of Sartorius and G-CON at**

Process Intensification Forum | September 17-19 | Göttingen, Germany

[Register now](#)

**Learn more about Sartorius ProcessGo®:**

<https://www.sartorius.com/en/pr/processgo>

**About Sartorius**

Sartorius is a leading international partner of 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 safely, rapidly, and economically. The company, based in Göttingen, Germany, 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 2023, the company generated sales revenue of around 3.4 billion euros. Currently, around 14,000 employees are working for customers around the globe.

Visit our [Newsroom](#) and follow Sartorius on [LinkedIn](#).

**About G-CON**

G-CON specializes in the design and construction of advanced cleanroom solutions for a wide range of industries. G-CON works closely with customers and A&Es to provide solutions that meet their specific needs, building PODular cleanrooms offsite, and providing modular and hybrid construction options. These offerings enable rapid deployment and easy configuration of cleanrooms while delivering the highest quality facilities, quickly, on time and on budget as well.

Connect with G-CON via [X](#) (formerly Twitter), [LinkedIn](#), and [Facebook](#)

**Sartorius Contact**

Lars Hanf

Head of Marketing Bioprocess Solutions

Sartorius

[lars.hanf@sartorius.com](mailto:lars.hanf@sartorius.com)

+49 551 308 4492

**G-CON Contact**

Brittany Berryman

Director of Marketing

G-CON

[bberryman@gconbio.com](mailto:bberryman@gconbio.com)

+1 979 431 0700