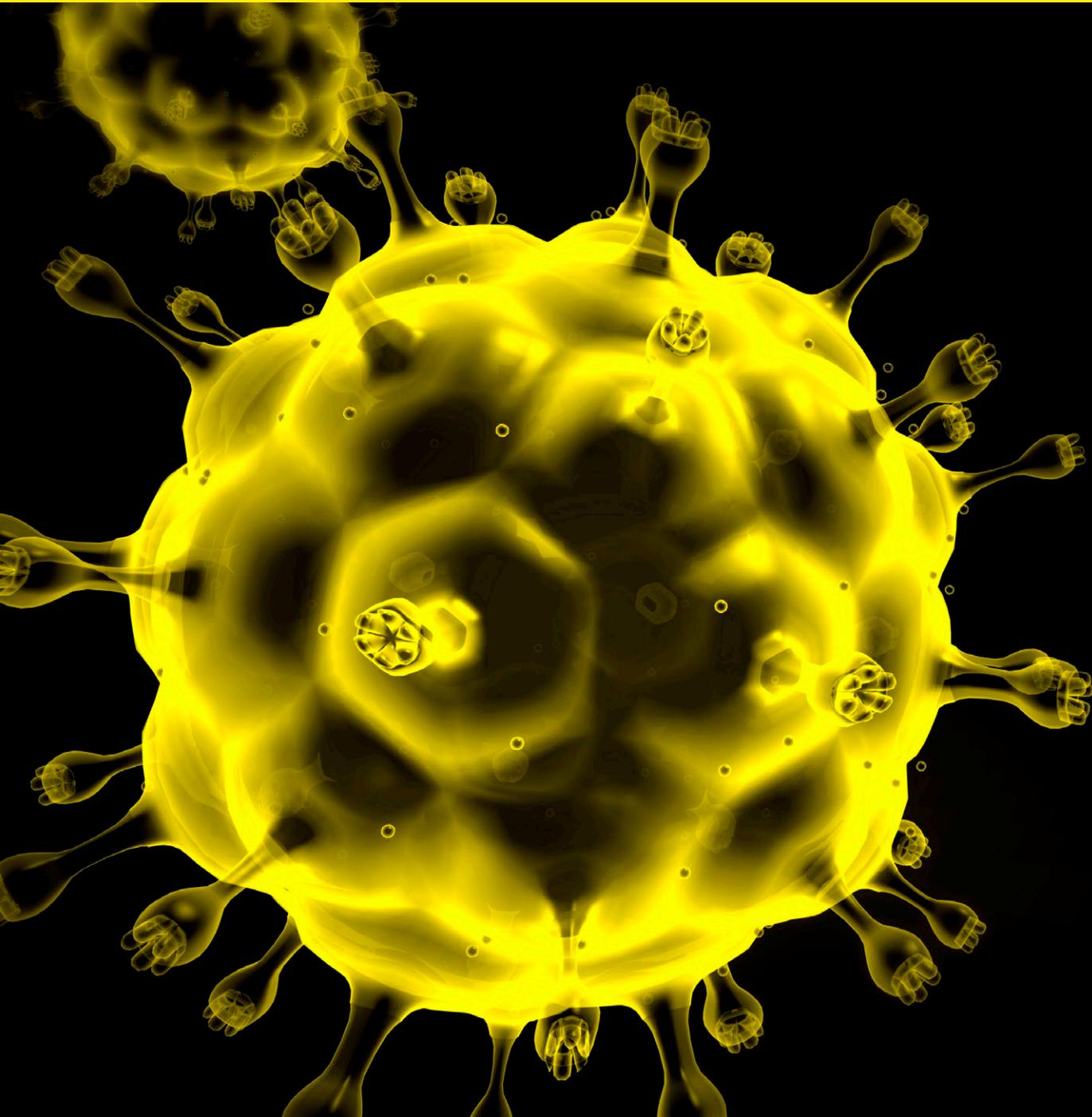


SARTORIUS

Emerging-Therapy-Services

Process Development Services for Novel Modalities



A Trusted Partner in Building Scalable and Efficient Processes for New Modalities From the Beginning

We are living in the century of biology; global medical knowledge continues to expand at a rapid pace. Disease breakthroughs continue, yet significant therapeutic needs remain unmet, such as treatments for cancers, rheumatic diseases, and Alzheimer's disease.

Sartorius is Part of the Solution

We help to ensure that new scientific discoveries can be translated quickly into effective patient care. Our technologies, scalable manufacturing platforms with advanced controls and analytics, coupled with our extensive partnership networks help achieve a common goal: better health for more people.

Simplifying Your Path to the Clinic

The cell and gene therapy market is growing strongly. While there have been some major breakthroughs, the industry faces significant challenges: developing scalable, robust processes, meeting regulatory demands, reducing costs, and speeding up time-to-market.

Sartorius has built an attractive portfolio of innovative technologies and services that streamline processes from research and development (R&D) to process development and manufacturing in the cell and gene therapy space. These solutions are designed to help biopharma customers deliver cell and gene therapies, as well as other new drug modalities, to patients more easily and cost-effectively. Whether it is process development or clinical phase manufacturing, our portfolio of proven solutions helps scientists achieve faster time-to-market while saving costs.

Leverage the experience and technical expertise to build efficient, scalable processes for your molecule.

- Get direct access to process scientists and the latest innovative technologies to support new and developing therapeutic modalities.
- De-risk your therapy with access to our extensive portfolio of products and services including solutions for cells, media, and pDNA.
- Gain knowledge of your molecule and process enabled by our expertise, inprocess analytics and hybrid (AI | ML) mechanist models.

State of the Art Facilities

- US-based facilities situation in Ann Arbor, MI & Marlborough, MA
- 15,000 sq ft to support multiple modalities
- EU Annex 1 Grade B & ISO Class 5 Cleanrooms
- Nine Process Development Suites



For further information, contact
tiffany.pogue@sartorius.com

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For more information, visit
sartorius.com