



SARTORIUS

Maximize Your Research with High-Quality Cell Culture Media, Reagents, and Supplements

Cell culture media is a nutrient-rich solution designed to support the growth, sustenance, and proliferation of cells in an artificial environment. It typically contains a blend of essential nutrients such as amino acids, vitamins, minerals, glucose, and growth factors, as well as buffering agents to maintain the pH at a level conducive to cell growth. The media can be tailored to the specific needs of different cell types, whether they are bacterial, plant, or animal cells.

The importance of cell culture media lies in its role in providing a controlled and replicable environment for cells outside of their natural context, such as within an organism or tissue. This is crucial for a variety of applications in life science research.

Sartorius offers a full portfolio of high-quality cell culture media, buffers, transfection reagents, growth factors, cytokines, and microtissue technologies for research applications so you can craft the perfect cell culture recipe for success in your lab.



Biological Research

It allows scientists to study cellular processes, genetics, and biochemistry in a detailed and controlled manner.



Drug Development

Cells grown in culture can be used for evaluating the efficacy and toxicity of new pharmaceutical compounds.



Vaccine Production

Cell cultures are used to grow viruses that are then used to produce vaccines.



Regenerative Medicine

Cultured cells are essential for tissue engineering and regenerative therapies.



Cancer Research

Cell culture enables the study of cancer cells in isolation, allowing for better understanding and treatment development.

Our Solutions

Research Grade Growth Factors and Cytokines

Sartorius offers a range of high-quality research use only (RUO) cytokines and growth factors which are produced using recombinant DNA technology and do not contain any animal-derived components or contaminants.

Our highly validated RUO Growth Factors and Cytokines use relevant assay models and workflows providing reliable and reproducible data and supply chain consistency. They are manufactured to the highest quality standards, ensuring high purity and efficacy, providing lot-to-lot consistency and low endotoxicity. Sartorius has implemented a quality management system certified for compliance with ISO9001.

- Animal-free and carrier protein-free
- Stringent validation using relevant assay models and workflows provides reliable and reproducible data
- Supply chain consistency enable seamless transition from research to production
- Maximum quality and safety reassurance due to state-of-the-art production, stringent control, and expert technical support



Classical Cell Culture Media and Reagents

Ensure quality cell culture with a wide range of classical media and reagents to support the growth and maintenance of a variety of cells and cell lines. From establishing and maintaining basic cell cultures to more advanced cell models, Sartorius offers a range of optimized ready-to-use products designed to meet your research needs. Each lot is manufactured under a strictly controlled process according to a Product Master Record to provide lot-to-lot consistency. Our excellence and expertise provide you with the safest, most dependable, and consistent cell culture media products.



Transfection Reagents for Life Science Research

Our extensive range of ready-to-use transfection reagents are suitable for research involving most mammalian primary cells and cell lines. These best-in-class transfection products allow scientists to study the function of a protein of interest, its role in a signaling pathway, perform genome editing using plasmid-based or RNA-based proven methods, and to conduct cost-effective RNA interference to transfect guide RNA and express Cas9 protein in mammalian cells.



Product name	SKU Number	Size
INTERFERin® siRNA/miRNA transfection	101000036	0.1 mL
	101000028	1 mL
	101000016	5x1 mL
PULSin® protein delivery	101000010	0.4 mL
jetMESSENGER® mRNA transfection	101000056	0.1 mL
	101000005	0.75 mL
jetOPTIMUS® DNA transfection	101000051	0.1 mL
	101000025	0.75 mL
	101000006	1.5 mL
jetOPTIMUS® buffer	201000001	60 mL
jetPEI® DNA transfection	101000053	1 mL
	101000020	4x1 mL
jetPRIME® DNA/siRNA (co-) transfection	101000027	0.1 mL
	101000015	0.75 mL
	101000046	1.5 mL
	101000001	5x1.5 mL
jetPRIME® buffer	201000003	60 mL

NexaGel® 3D Cell Culture Matrices

*NexaGel® hydrogel is a ready-to-use, versatile synthetic matrix system that closely replicates the human microenvironment, making it ideal for research involving a wide range of cell types. NexaGel® hydrogels feature optimized multi-functional ligands and concentration formulations, suitable for various applications, including 3D cell models, stem cell spheroids, and organoids.



*NexaGel® is a registered trademark of Sartorius Bioanalytical Instruments, Inc. For details on the registrations please refer to our website www.sartorius.com/en/patents-and-trademarks.

Microtissue Technologies, Primary Cells, Cultureware, and Testing Services

Explore ready-to-use 3D human-derived microtissue models, human primary cells, high-quality glass bottom cell culture dishes and plates for superior imaging, as well as toxicology testing services for industries such as cosmetics, pharmaceutical, chemical, and other regulated industries.

🌐 Discover Microtissue Technologies



Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 650 322 1360
Toll-free +1 800 635 2906

🌐 For further information, visit
www.sartorius.com