



Comprehensive
Media Solutions
for Cell and Gene
Therapies

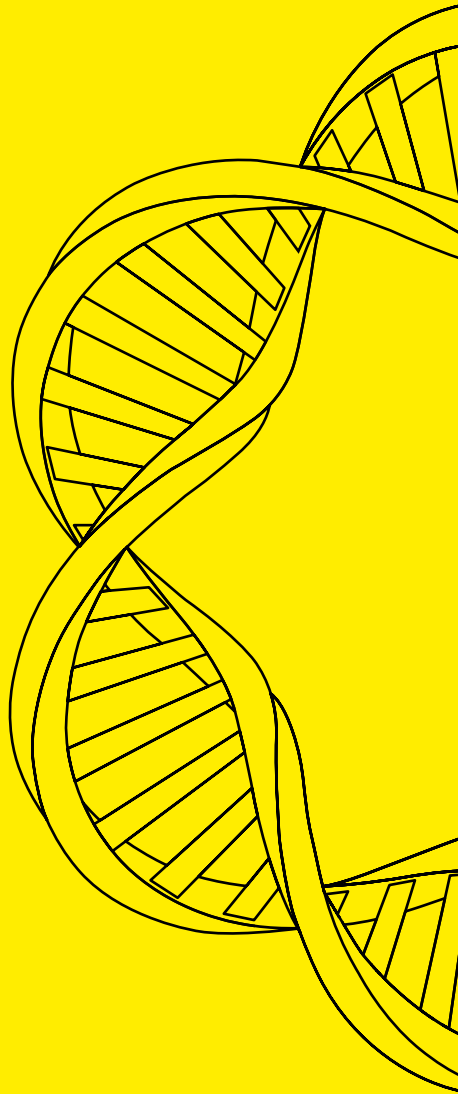
Simplifying Progress

SARTORIUS

Sartorius as Your Media Solutions Partner

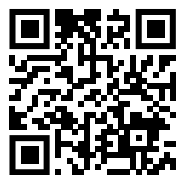
Maximize cell quality attributes,
ensure reproducible results, and reduce
your development timelines.

We offer high quality cell culture media
and reagents, critical raw materials,
biopreservation solutions, as well as a
variety of supporting services to cover
all major cell types in cell and gene
therapies. Backed by expert support
and a global manufacturing network,
we provide what's vital to your cells
and your success – from screening to
commercial production.



Request a Sample

sartorius.com/request-sample-cgt







Stem Cell Media, Supplements, and Reagents

Human Induced Pluripotent Stem Cells | Embryonic Stem Cells (hiPSCs/hESC)s

Serum-free, xeno-free media portfolio for iPSC and ESC culture. Manufactured ISO 13485 in compliance with applicable cGMP guidelines, DMF available.

Product	Format	Features
NutriStem® hPSC XF Medium w/ or w/o Phenol Red	Bottle (500 mL), bag (1 L, 1.5 L, 2 L, 2.8 L, 10 L)	Specially formulated for optimal growth and expansion of pluripotent stem cells
NutriStem® hPSC XF Medium w/o TGF & FGF	Bottle (500 mL)	Growth factor-free formulation for the growth and expansion of PSCs
NutriStem® hPSC XF Medium with +20 ng/mL bFGF	Bottle (500 mL)	Modified formulation containing +20 ng/mL bFGF ideal for adaptation of PSC cultures from serum-containing or other commercial medium

See Page 7 to 10 for additional reagents that support hiPSCs/hESC)s

Human Mesenchymal Stromal Cells (MSCs)

Serum-free, xeno-free culture media and substrate portfolio for MSC isolation, expansion and differentiation. Manufactured ISO 13485 in compliance with applicable cGMP guidelines, DMF available for NutriStem.

Product	Format	Features
MSC NutriStem® XF Basal Medium w/ or w/o Phenol Red	Bottle (100 mL, 500 mL), bag (5 L, 10 L, 20 L)	Basal medium designed for optimal growth and expansion of MSCs derived from multiple sources. Intended for use with MSC NutriStem XF Supplement
MSC NutriStem® XF Supplement	Bottle (30 mL, 60 mL, 120 mL), vial (0.6 mL, 3 mL, 9 mL)	Supplement formulated for optimal growth and expansion of MSCs derived from multiple sources. Intended for use with MSC NutriStem XF Basal Medium
MSCgo™ Chondrogenic Differentiation Medium	Bottle (100 mL)	Basal medium optimized for directed differentiation of MSCs to chondrocytes when used with MSCgo Chondrogenic Differentiation Supplement
MSCgo™ Chondrogenic Differentiation Supplement Mix	Vial (10 mL)	Supplement optimized for directed differentiation of MSCs to chondrocytes when used with MSCgo Chondrogenic Differentiation Basal Medium
MSCgo™ Adipogenic Differentiation Medium	Bottle (100 mL)	Basal medium for directed differentiation of MSCs to adipocytes when used with MSCgo Adipogenic Differentiation Supplement Mix 1 & 2
MSCgo™ Adipogenic SF XF Supplement Mix 1	Vial (0.1 mL)	Supplement optimized for directed differentiation of MSCs to adipocytes when used with MSCgo Adipogenic Differentiation Medium & Supplement Mix 2
MSCgo™ Adipogenic SF XF Supplement Mix 2	Vial (1.5 mL)	Supplement optimized for directed differentiation of MSCs to adipocytes when used with MSCgo Adipogenic Differentiation Medium & Supplement Mix 1
MSCgo™ Osteogenic Differentiation Medium	Bottle (100 mL)	Complete medium optimized for directed differentiation of MSCs to osteoblasts
MSCgo™ Rapid Osteogenic Differentiation Medium	Bottle (100 mL)	Complete medium optimized for faster directed differentiation of MSCs to osteoblasts
NutriCoat® Solution	Vial (1.5 mL)	Defined substrate based on clinical-grade human Fibrinogen, designed for the attachment of MSCs in serum-free and xeno-free culture systems

See Page 7 to 10 for additional reagents that support MSCs

Human Hematopoietic Stem Cells (HSCs)

Serum-free, xeno-free culture medium for expansion of HSCs. Manufactured in accordance with applicable GMP guidelines. DMF and PDMA available.

Product	Format	Features
CellGenix® GMP SCGM w/ or w/o Phenol Red	Bottle (500 mL), bag (500 mL)	Optimized for serum-free expansion of low numbers of isolated human HSC and progenitor cells (CD34+ cells)

See Page 7 to 10 for additional reagents that support HSCs

Immune Cell Media

T Cells

Serum-free, xeno-free T cell and immunotherapy culture media portfolio optimized for culturing PBMC, CAR-T, TIL, Treg. Manufactured ISO 13485 in compliance with applicable cGMP guidelines.

Product	Format	Features
4Cell® Nutri-T GMP Medium w/ or w/o Phenol Red	Bottle (1 L), bag (1 L, 5 L, 10 L)	Formulated for the culture and expansion of lymphocytes for cell-based immunotherapy, tested and developed on both healthy and non-healthy patient derived donor cells
CellGenix® GMP TCM	Bottle (500 mL), bag (1 L)	Formulated for serum-free expansion and cultivation of human T cells, highly functional CAR-T cells, and Tregs, in static or simple stirred bioreactors. DMF available.

See Page 7 to 10 for additional reagents that support T cells

Natural Killer Cells (NK Cells)

Serum-free, xeno-free culture medium for expansion and cultivation of NK cells. Manufactured in accordance with applicable GMP guidelines. DMF and PDMA available for CellGenix GMP SCGM.

Product	Format	Features
CellGenix® GMP SCGM w/ or w/o Phenol Red	Bottle (500 mL), bag (500 mL)	Optimized for serum-free expansion of low numbers of isolated human HSC and progenitor cells (CD34+ cells). Highly efficient for the expansion and cultivation of human NK cells. DMF and PDMA available.
4Cell Nutri-T GMP Medium w/ or w/o Phenol Red	Bottle (1 L), bag (1 L, 5 L, 10 L)	Formulated for the culture and expansion of lymphocytes (including NK cells) for cell-based immunotherapy, tested and developed on both healthy and non-healthy patient derived donor cells

See Page 7 to 10 for additional reagents that support natural killer cells

Dendritic Cells (DCs)

Serum-free, xeno-free culture medium for the generation of DCs. Manufactured in accordance with applicable GMP guidelines. DMF available.

Product	Format	Features
CellGenix® GMP DC w/ or w/o Phenol Red	Bottle (500 mL), bag (500 mL)	Optimized to support the serum-free differentiation from CD14+ monocytes to immature DCs or macrophages, and further from immature to mature DCs.

See Page 7 to 10 for additional reagents that support dendritic cells



Recombinant Human Albumin

Ready-to-use, animal-component free, serum-free, chemically-defined, manufactured to ICH Q7 cGMP standards, extensive DMF available.

Product	Format	Applications
The Recombumin® Portfolio	Vial (50 mL), bioprocess container* (1 L)	Supports iPSCs ESCs, MSCs, HSCs, T cells, NK cells, DC; Applications include upstream bioprocessing, cell culture supplement, downstream processing, cryopreservation and final formulation excipient

*Selected products

Biopreservation Solutions

Cryopreservation and Cold Storage Solutions

Chemically defined, animal component-free cryopreservation and cold storage solutions designed and validated for the preservation of sensitive cells such as stem cells and immune cells. Manufactured ISO 13485 in compliance with applicable cGMP guidelines, DMFs available.

Product	Format	Features
NutriFreez® D10 Cryopreservation Medium, w/ or w/o Phenol Red	Bottle (10 mL, 20 mL, 50 mL, 100 mL, 500 mL)	Ready-to-use cryopreservation medium containing 10% DMSO
NutriFreez® D5 Cryopreservation Solution, w/o Phenol Red	Bottle (10 mL, 100 mL, 500 mL), bag (100 mL)	Ready-to-use, salt-based cryopreservation solution containing 5% DMSO
NutriStor® Cold Storage Solution	Bottle (10 mL, 100 mL, 500 mL), bag (250 mL, 500 mL)	Designed for short-term storage and shipment of highly sensitive cells at 2 - 8 °C for clinical applications



Other Media and Reagents

Endothelial Cell Media

Serum-free, xeno-free culture media for endothelial cell expansion. Manufactured ISO 13485 in compliance with applicable cGMP guidelines.

Product	Format	Features
EndoGo™ XF Basal Medium	Bottle (500 mL)	Basal medium designed for long-term expansion of macro- and microvascular endothelial cells. Intended for use with EndoGo XF Supplement
EndoGo™ XF Supplement	Vial (2.5 mL)	Supplement designed for long-term expansion of macro- and microvascular endothelial cells. Intended for use with EndoGo XF Basal Medium

Other Supplements

Product	Format	Features
Bio-Pure Human Serum Albumin 10% Solution	Bottle (10 mL, 100 mL)	Specially treated Human Serum Albumin, optimized for ES and iPS cell culture
Human Serum Albumin 20% Solution	Bottle (50 mL, 500 mL)	Human Serum Albumin for ES and iPS cell culture
PLTGold® Human Platelet Lysate (Clinical Grade)	Bottle (27 mL, 100 mL, 500 mL, 1 L)	Alternative to FBS for MSC, HSC, PSC, ESC, EC and other primary cell culture. Gamma irradiated, US origin
PLTGold® Human Platelet Lysate (GI Clinical Grade)	Bottle (27 mL, 100 mL, 500 mL, 1 L)	Alternative to FBS for MSC, HSC, PSC, ESC, EC and other primary cell culture. Clinical-grade, US origin
PLTGold® Human Platelet Lysate (Research Grade)	Bottle (27 mL, 100 mL, 500 mL, 1 L)	Alternative to FBS for MSC, HSC, PSC, ESC, EC and other primary cell culture. Research-grade, US origin

Dissociation Solutions

Product	Format	Features
Recombinant Trypsin Solution	Bottle (20 mL, 100 mL)	Animal component-free cell dissociation solution
Recombinant Trypsin EDTA Solution	Bottle (20 mL, 100 mL, 500 mL)	Animal component-free cell dissociation solution with accelerated enzymatic action



Growth Factors and Cytokines

Animal- and human-component free, preclinical and GMP (if applicable), DMF available.

Product	Format	Applications
CellGenix® rh Activin A* (Preclinical and GMP Grade)	Vial (50 µg)	Cell culture for iPSCs ESCs
CellGenix® rh EGF (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for iPSCs ESCs, and MSCs,
CellGenix® rh FGF-2 (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for iPSCs ESCs, and MSCs
CellGenix® rh Flt-3L (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for HSCs, and NK
CellGenix® rh GM-CSF (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for DC
CellGenix® rh IFN-γ* Preclinical grade only	Vial (50 µg)	Cell culture for DC
CellGenix® rh IL-10 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for DC
CellGenix® rh IL-15 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for T cells, and NK cells
CellGenix® rh IL-1β (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for iPSCs ESCs and DC
CellGenix® rh IL-2 (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for T cells, and NK cells
CellGenix® rh IL-21 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for T cells, and NK cells
CellGenix® rh IL-3 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for HSCs
CellGenix® rh IL-4 (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for DC
CellGenix® rh IL-6 (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for HSCs, NK cells, and DC
CellGenix® rh IL-7 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for T cells, and NK cells
CellGenix® rh OSM* Preclinical grade only	Vial (50 µg)	Cell culture for iPSCs ESCs
CellGenix® rh PDGF-BB* (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for MSCs
CellGenix® rh SCF (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for iPSCs ESCs, HSCs, and NK cells
CellGenix® rh TGF-β1 (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for iPSCs ESCs, MSCs, and T cells
CellGenix® rh TNF-α (Preclinical and GMP grade)	Vial (50 µg)	Cell culture for DC
CellGenix® rh TPO (Preclinical and GMP grade)	Vial (50 µg, 1 mg**)	Cell culture for iPSCs ESCs, HSCs, and NK cells

* DMF currently not available

** 1 mg format only available for GMP-grade cytokines



HEK293 Media Portfolio

Chemically-defined, animal-component free, serum-free, protein-free, and hydrolysate-free media and feed portfolio for HEK293 suspension culture. These media are designed for viral vector production and support high and consistent titers in both clinical and commercial applications.

Product	Format	Features
HEK GM CD Medium	Bottle (1 L), bag (10 L*, 20 L, 50 L, 100 L, 200 L, 500 L), powder (10 L*, 100 L, 500 L)	With growth factor, a robust and effective base medium for infection and stable expression
HEK TF CD Medium	Bottle (1 L), bag (10 L*, 20 L, 50 L, 100 L, 200 L, 500 L), powder (10 L*, 100 L, 500 L)	With growth factor, a well-established transfection medium for growth, transfection, and infection
HEK FS_2 Feed	Bottle (1 L), bag (5 L, 10 L, 20 L), powder (5 L, 10 L, 50 L)	Boost viral production for gene therapies and vaccines. This feed can be used with the 4 HEK media portfolio.
HEK ViP NB CD Medium	Bottle (1 L), bag (10 L*, 20 L, 50 L, 100 L, 200 L, 500 L), powder (10 L*, 100 L, 500 L)	Without growth factor, the leanest formulation with high performance, for growth, transfection, and infection
HEK ViP NX CD Medium	Bottle (1 L), bag (10 L*, 20 L, 50 L, 100 L, 200 L, 500 L), powder (10 L*, 100 L, 500 L)	Without growth factor, the highest nutrient level content for demanding cells, for growth, transfection, and infection. Little or no adaptation required from other serum-free media

*in stock



Media Optimization Services For Viral Vector Production

We are equipped to meet your demand in cell culture products and services regardless of the current workflow stage. We offer Media Customization Services for our catalog products (add or remove components, selection of formats, customized packaging options, QC testing selection).

We leverage our resources to design robust cell culture media tailored to your specific cell line requirements for consistent performance. From small scale non-GMP manufacturing in our Media Rapid Prototyping line to commercial batches in our animal origin-free plant, we partner with you to support your entire cell culture journey.

- Optimization & Development Services for Feed and Media
- Cell Line Adaptation Service
- Benchmarking & Material Testing
- Tech Batches & Feasibility Studies
- Media Rapid Prototyping
- Large scale manufacturing service

Media and Spent Media Analysis Service

Spent media analytics is a valuable tool that enables the tracking of changes in medium composition. The information gathered reveals insights into the metabolic processes of the cell population and how the media influences process and product characteristics. Services include:

- ELISA-based analysis for AAV (2, 5, 8, 9)
- Custom method establishment
- Custom analytical measurements
- Targeted and untargeted analytics
- Amino Acid Analysis
- Dipeptide Analysis by UHPLC-MS/MS
- Analysis of Total Amino Acid Content
- Water-Soluble Vitamins by UHPLC-MS/MS
- Sucrose Analysis
- Ascorbate (Vitamin C) Analysis
- Glucose and Sucrose Analysis
- Glutathione (oxidized & reduced) Test
- Phosphate Analysis by Photometric assay
- Glucose Analysis
- Glucose and Lactate Analysis
- Ammonium (Concentration) Analysis
- Pyridoxal Phosphate Measurement by UHPLC
- Culture Media Analysis Bundle
- High Abundant Cation Analysis by ICP-MS
- Organic Acids Analysis by UHPLC-MS
- Peptide Size Determination by SEC
- Spent Media – Trace Element Bundle
- Polyamine Analysis by UHPLC-MS/MS
- Element Analysis Bundle (5, 10, 15, 20, or 25 elements)

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 further information, visit**
[sartorius.com](https://www.sartorius.com)