



High Performance Cell Culture Media, and Reagents, and Supplements for Life Science Research



New

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



Transfection Reagents for Life Science Research

- Extensive range of ready to use transfection reagents suitable for research involving most mammalian cell lines
- 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
- Conduct cost effective RNA interference to transfect guide RNA and express Cas9 protein in mammalian cells.



Research Grade Growth Factors and Cytokines

- Fully animal derived and component free
- Stringent validation guarantees reliable and reproducible data
- Supply chain consistency enable seamless transition from research to production



Classical Cell Culture Media and Reagents

- Wide range of high quality media and reagents to support the growth and maintenance of a variety of cells and cell lines
- Strictly controlled manufacturing process provides lot to lot consistency



NexaGel® High Performance Hydrogel

- Xeno-free: 100% animal origin-free
- Ready-to-use formulation: No activation or adjustments required
- Easy and fast cell harvesting: Recover cells in just 20 minutes



Elevate your research: Try a sample of our research grade Cytokines or Transfection Reagents!



Research Grade Growth Factors
and Cytokines



Scan to submit
your sample
request:



Transfection Reagents for Life
Science Research



Scan to submit
your sample
request:



For testing of NexaGel® and Classical
media please contact your local
Sartorius customer service or sales
colleague.

For more information:

Additional products supporting
advanced research in pre-clinical
and GMP areas.



Germany

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

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

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