



Comprehensive
Media Solutions for
Your Protein-Based
Therapies

Simplifying Progress

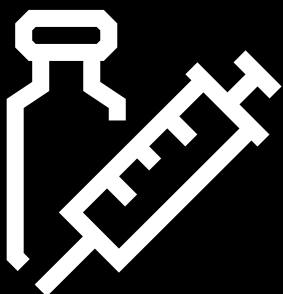
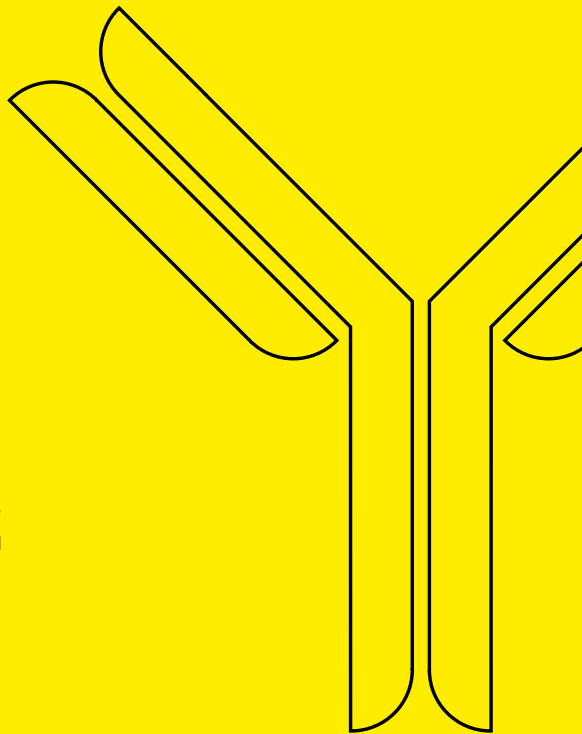
SARTORIUS

Sartorius as Your Media Solutions Partner

Chinese hamster ovary (CHO) cells are ideal for therapeutic protein production in biopharmaceutical research and development.

Our comprehensive offering of several CHO cell culture media and companion feeds, combined with our advanced cell line development and supporting media services, enables simplified clone selection and optimized titer production.

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

[www.sartorius.com/en/pr/
request-cell-culture-media-sample](http://www.sartorius.com/en/pr/request-cell-culture-media-sample)





4Cell® CHO Platform

The 4Cell® CHO Platform bundles services and media in a GMP manufacturing-ready solution for drug developers looking for stable and ready-to-scale clones.

Cell Line Development Service

The platform comprises an advanced CLD service using Sartorius' high-throughput CellCelector™ technology for Single Cell Cloning that can deliver high-performing clones in 9 weeks with demonstrated performance and titers up to 10 g/L in the optimized 4Cell® SmartCHO media.

Cell Culture Media

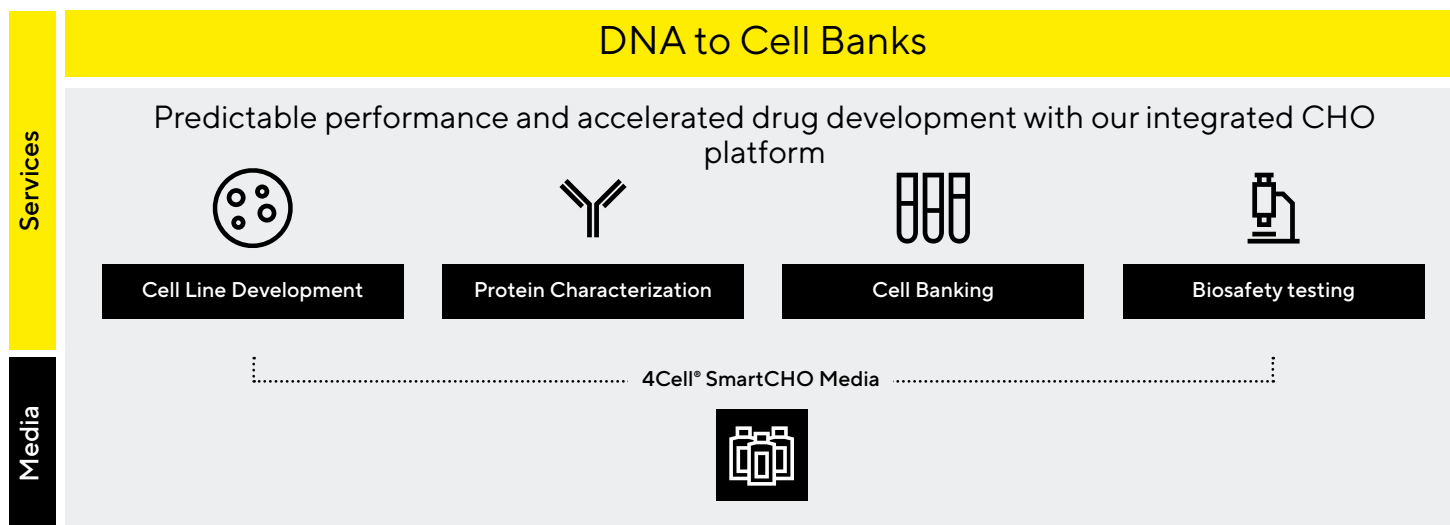
The 4Cell® SmartCHO Media was specially developed for our cell line development service and supports the production of RCB and MCB. It is a high-quality, high-performance cell culture formulation that is chemically defined and free of animal components to maximize product titer in our CHO cell line. 4Cell® SmartCHOpe Media, our perfusion medium, is compatible with our CLD service and supports clone selection in a perfusion mimic protocol.

Protein Characterization Service

Sartorius' protein characterization services are scaled to each phase of drug development. These services include the development, qualification, and validation of product characterization assays, such as cell-based potency assays, binding assays, and physicochemical tests.

Cell Bank Manufacturing And Biosafety Testing

The Sartorius cell banking platform is designed as a closed, sterile, single-use system with automated filling into vials to generate large and consistent banks. Plus, Sartorius has extensive experience in characterizing cell banks used in the production of protein-based therapies.



CHO Media and Feeds

Sartorius offers a wide range of media and feeds to provide drug developers with the best formulation for their stable clones, enabling a successful transition to the clinic with high-performance cell growth and productivity.

Chemically defined, animal ingredient-free, protein-free, serum-free, hydrolysate-free, and GMP-compliant media to support the expansion of the most common CHO cell lines in all applications.

CHO Transfection Media

Product	Format	Features
CHO TF	Bottle (1 L), powder (10 L, 50 L)	Ideal medium for working with transient CHO cells

CHO Batch and Fed-Batch Media

Product	Format	Features
4Cell® SmartCHO SAM	Bottle (1 L), bag (1 L, 5 L, 20 L), powder (10 L, 25 L, 250 L)	Allows the cultivation or media adaptation of CHO cells
4Cell® SmartCHO PM	Bottle (1 L), bag (1 L, 5 L, 20 L), powder (9 L, 20 L, 200 L)	Non-selective production media for batch or fed-batch culture. It can be used for effective cultivation of any CHO cell lines, and helps produce high yield of the active protein
4Cell® SmartCHO FMA	Bottle (1 L), bag (1 L, 5 L, 20 L), powder (3 L, 30 L)	Ideal companion feeds designed to be used together with 4Cell® SmartCHO Production Media
4Cell® SmartCHO FMB	Bottle (125 mL, 500 mL), bag (1 L, 5 L), powder (4 L, 45 L)	Ideal companion feeds designed to be used together with 4Cell® SmartCHO Production Media
CHOlean	Bottle (1 L), powder (10 L, 50 L, 100 L)	Lean and robust growth media for a variety of applications, lean formulation with minimal components ideal for media optimization activities
TC-42	Bottle (1 L), powder (10 L, 50 L)	Workhorse formulation especially designed for CHO-DHFR cell lines
TCX10D	Bottle (1 L), powder (10 L, 50 L, 100 L, 200 L, 500 L)	For long-term high-performance growth with CHO-GS cell lines
TCX6D	Bottle (1 L), powder (10 L, 50 L, 100L, 200 L, 500 L)	One high-performing media for any CHO clones with growth factor LONG® R3IGF-I
TCX7D Feed Supplement	Bottle (1 L), powder (2 L, 5 L, 10 L, 25 L, 50 L, 100 L)	With 60 g/L D-glucose, especially for CHO GS cell lines
Basic Feed	Bottle (1 L), bag (10 L, 20 L, 50 L), powder (5 L, 10 L, 50 L)	With 20 g/L D-glucose, also suitable for hybridoma



CHO Perfusion Media

Product	Format	Features
4Cell® SmartCHO SAM	Powder (10 L, 25 L, 250 L)	Allows the cultivation or media adaptation of CHO cells
4Cell® SmartCHO PM	Powder (20 L, 50 L, 200 L, 900 L)	Non-selective production media for perfusion culture
4Cell® SmartCHO FMA	Powder (3 L, 30 L, 120 L)	Ideal companion feeds designed to be used together with 4Cell® SmartCHOpe Production Media
4Cell® SmartCHO FMB	Powder (2 L, 4.5 L, 45 L, 180 L)	Ideal companion feeds designed to be used together with 4Cell® SmartCHOpe Production Media



Media Optimization Services

Supporting customers in optimizing their media and processes to prepare for scale-up, decreasing their time-to-market while maintaining high titer and product quality

- Cell Line Adaptation Service
- Media and Process Optimization Service
- Benchmarking and Material Testing
- Tech Batches and Feasibility Studies
- Spent Media Analysis

Custom Media Development And Manufacturing Services

Targeted, rapid, and efficient services for the development of chemically defined, animal component-free, custom cell culture media and feed, available as liquid or powder formulations.



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**
www.sartorius.com