

4Cell® XtraCHO Media System

Large Scale: Media System
for Protein Production
Using CHO Systems

Benefits

Use of production media and feeding supplements helps satisfying cell culture requirement using different approaches conditions. Recommended protocols facilitate easy and efficient identification of a feeding strategy.

- **Production Efficacy:** Ideal from bench scale to manufacturing
- **Compliance with Regulatory Guidelines:** Eliminate process variability and contamination risks
- **Consistency and Traceability:** Fully chemically defined medium
- **Reduce Process Complexity:** Simplified purification

Product Information

The Starting Kit of the 4Cell® XtraCHO Media System includes all the necessary media components from adapting the cells to fed-batch production at a cultivation volume of 3L. For larger volumes all individual system components are available in different formats, in liquid and powder.

Stock & Adaptation Medium: It allows the cultivation or media adaptation of cells. Lacking certain metabolites, the cultivation media exerts a strong selection pressure on cells and shortens the time consuming adaptation to fed batch culture.



Production Medium: It is a non-selective production medium for batch or for batch culture. It can be used for effective cultivation of recombinant CHO DG44 cells, this media helps produce high yield of the active protein.

Feed Medium A and Feed Medium B: Both feed additives are especially matched to the production medium. Combining those increases fed-batch productivity 3- to 5-fold compared to batch culture conditions.

Applications

Protein and mAb production in DHFR deficient CHO DG44 cell lines and other CHO cell lines.

Features of the Standard Product



Technical Data

Specifications

Media Type	Chemically-defined medium: does not contain serum or undefined components
Easy-to-use	Adaptation, production and feeds all designed to work together
Cell Line	Developed for CHO DG44 cells and works also with other CHO-Systems
Storage Condition	2–8°C, protect from light
Shelf Life	12 months for liquid, 3 years for powder from date of manufacture

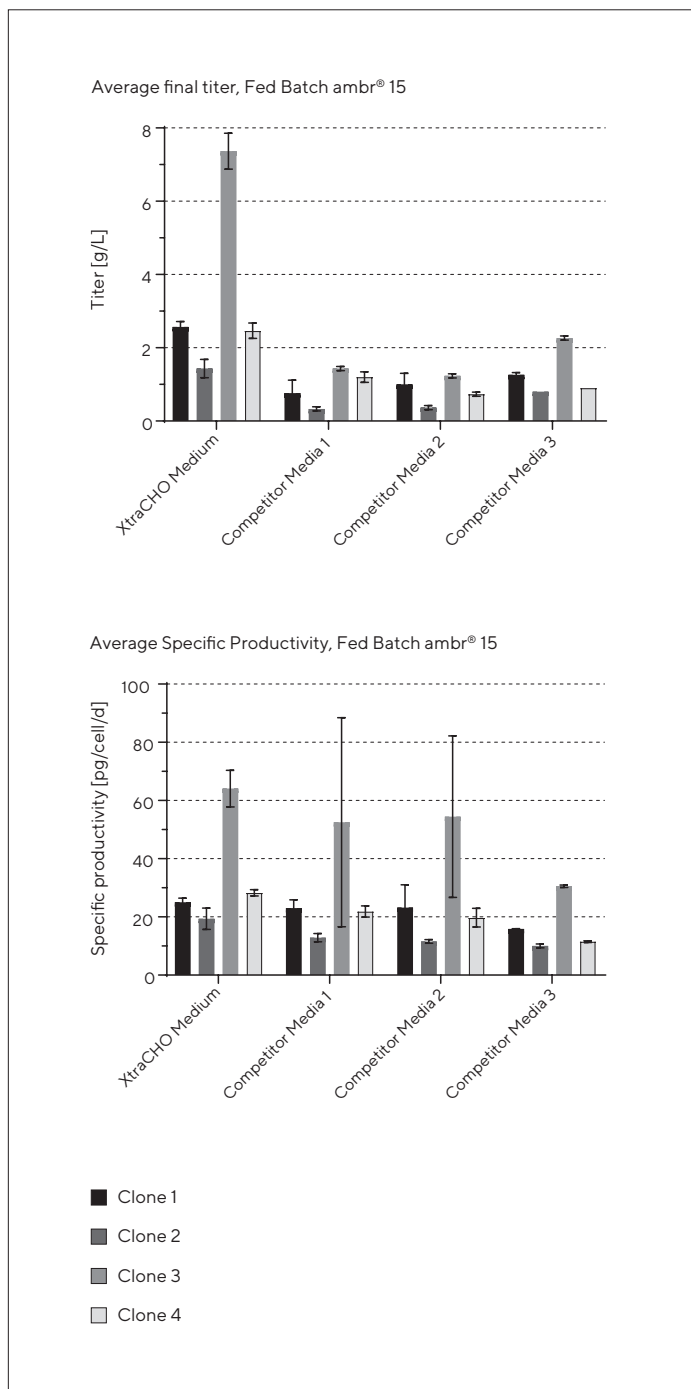
Features of the Standard Product

- CD** **Chemically Defined:** The exact concentration and size of every component is known
- NAO** **Non-Animal Origin:** the formulation is entirely made from non-animal | -human origin components
- PF** **Protein-free:** the formulation does not contain any protein
- FFM** For research or for further manufacturing use
- Product available in powder format
- Product available in liquid format

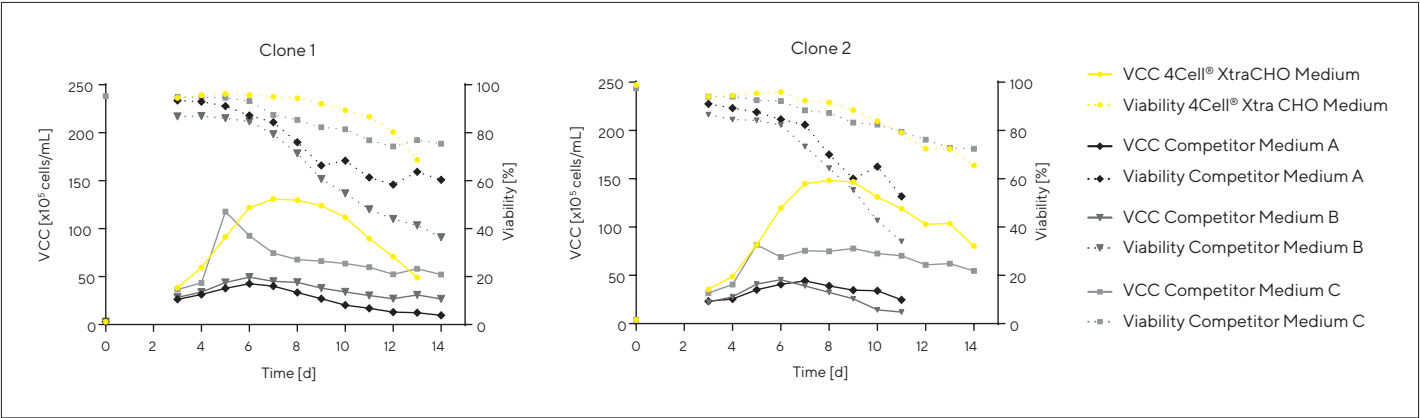
Performance

High product titer processes by maintaining high cell viability and high cell density throughout 14 days cultivation

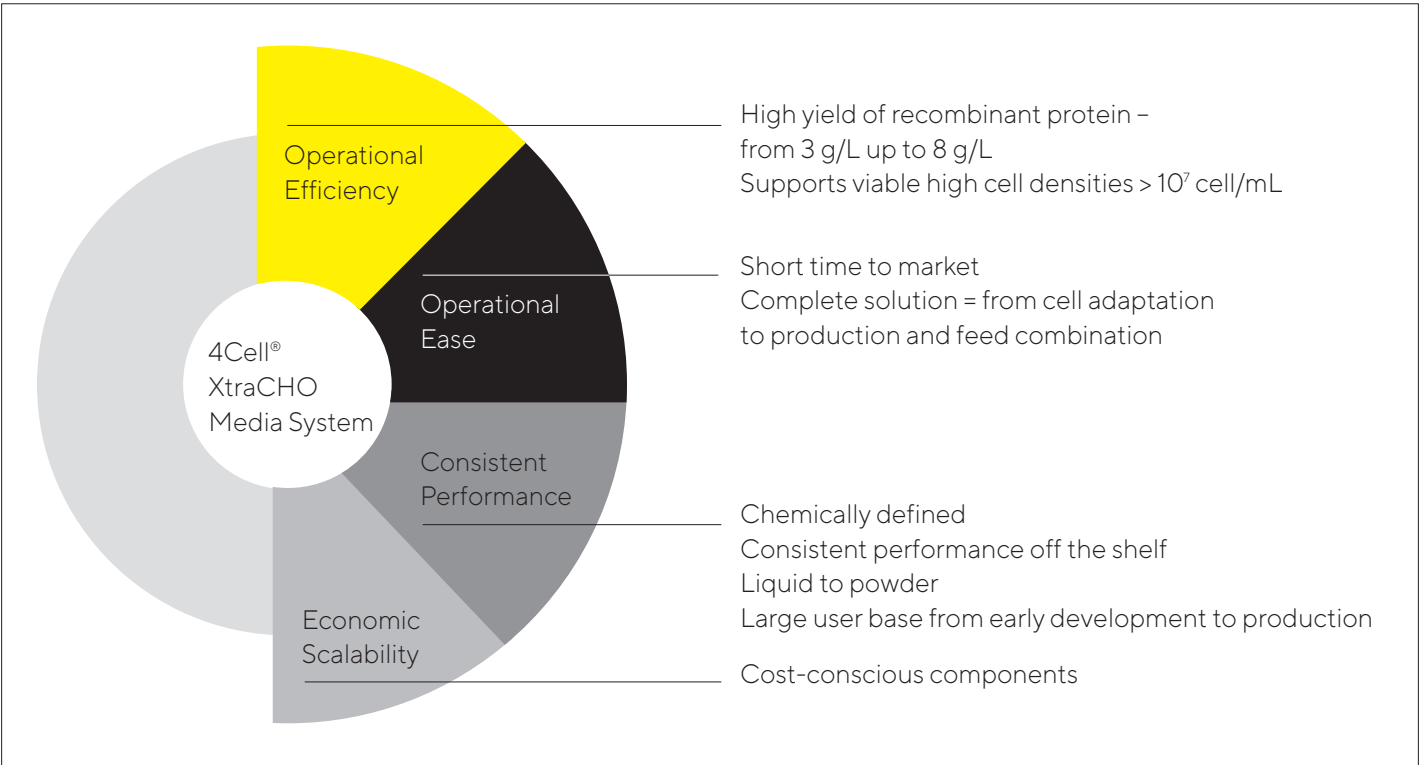
ambr® 15 Titer and specific productivity of 4 CHO DG44 Clones in XtraCHO Media System and 3 media of competitors in fed-batch mode



ambr® 250 Viable cell Concentration and Viability of 2 CHO DG44 Clones in 4Cell® XtraCHO Media System and 3 media of competitors in fed-batch mode



Your Benefits at a Glance



Ordering Information*

Description	Form	Package	Volume Size	Order Code
4Cell® XtraCHO Media System Starting Kit	Liquid	Bottle	125 – 1000 mL Kit (up to 3 L fed-batch)	CFP3FA0200
4Cell® XtraCHO Stock & Adaptation Medium (SAM)	Liquid	Bottle	1000 mL 2 units	CFP3FA1201
	Liquid	Bottle	1000 mL 6 units	CFP3FA1206
	Liquid	Bag	5 L 2 units	CFP3FA1101
	Powder	Bucket	3 L [60g] 1 unit	CQP3FA1114
	Powder	Bucket	15 L [300g] 1 unit	CQP3FA1113
	Powder	Bucket	25L [510g] 1 unit	CQP3FA1135
	Powder	Bucket	250L [5050g] 1 unit	CQP3FA1136
4Cell® XtraCHO Production Medium (PM)	Liquid	Bottle	1000 mL 2 units	CFP3FA2202
	Liquid	Bottle	1000 mL 6 units	CFP3FA2207
	Liquid	Bag	5 L 2 units	CFP3FA2104
	Liquid	Bag	20 L 1 unit	CFP3FA2105
	Powder	Bucket	6 L [135g] 1 unit	CQP3FA2115
	Powder	Bucket	30 L [670g] 1 unit	CQP3FA2116
	Powder	Bucket	200 L [4500g] 1 unit	CQP3FA2138
4Cell® XtraCHO Feed Medium A (FMA)	Liquid	Bottle	1000 mL 2 units	CFP3FA3203
	Liquid	Bottle	1000 mL 6 units	CFP3FA3208
	Liquid	Bag	5 L 2 units	CFP3FA3110
	Powder	Bucket	3 L [465g] 1 unit	CQP3FA3120
	Powder	Bucket	15 L [2310g] 1 unit	CQP3FA3121
	Powder	Bucket	30 L [4700g] 1 unit	CQP3FA3140
4Cell® XtraCHO Feed Medium B (FMB)	Liquid	Bottle	125 mL 2 units	CFP3FA4204
	Liquid	Bottle	500 mL 2 units	CFP3FA4205
	Liquid	Bottle	500 mL 6 units	CFP3FA4210
	Powder	Bucket	1 L [95g] 1 unit	CQP3FA4119
	Powder	Bucket	2 L [190g] 1 unit	CQP3FA4118
	Powder	Bucket	5L [480g] 1 unit	CQP3FA4142

* Other product formats available. For more information please contact:
 Europe | Asia: CellCultureMedia.EU@sartorius.com
 Americas: CellCultureMedia.NA@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 further contacts, visit
www.sartorius.com