

4Cell® BHK-21 CD Medium

Chemically Defined (CD) medium for BHK-21 cells for maximum cell growth and virus production

Benefits

Developed to maximize BHK-21 cell growth for virus and vaccine manufacturing, the chemically defined medium provides equivalent or higher growth of BHK-21 cells when compared to serum-supplemented conditions and offers:

- **Production Efficacy:** Ideal from bench scale to manufacturing
- **Serum-Free and Protein-Free:** Simplify your downstream purification and filtration processes
- **Non-Animal Origin and Chemically Defined:** Eliminate process variability and contamination risks
- **Antibiotic-free:** Eliminate risk of antibiotic traces



Product Information

The 4Cell® BHK-21 CD Medium is a chemically defined, serum-free, protein-free, animal component-free, hydrolysate-free medium designed to maximize the growth of BHK-21 cells in suspension culture. The 4Cell® BHK-21 CD Medium supports cell growth up to 7×10^6 cells/mL with greater than 90% viability. BHK-21 cells have been carried for more than 90 generations (30 passages) in 4Cell® BHK-21 CD Medium with no loss of viability.

This formulation is easy-to-use and is suitable for suspension cultures. It supports the superior production of both virus and recombinant proteins and requires only minimal adaptation from other serum-free media.

Applications

4Cell® BHK-21 CD Medium is a chemically-defined medium dedicated to cell growth, maximum virus propagation and protein expression. The medium is ideally suitable for production of a wide range of viruses using BHK-21 suspension cells:

- Foot & mouth
- Rabies for veterinary use
- Blue tongue
- Aujeszky's virus
- New castle and Polio vaccines

Features of the Standard Product



Technical Data

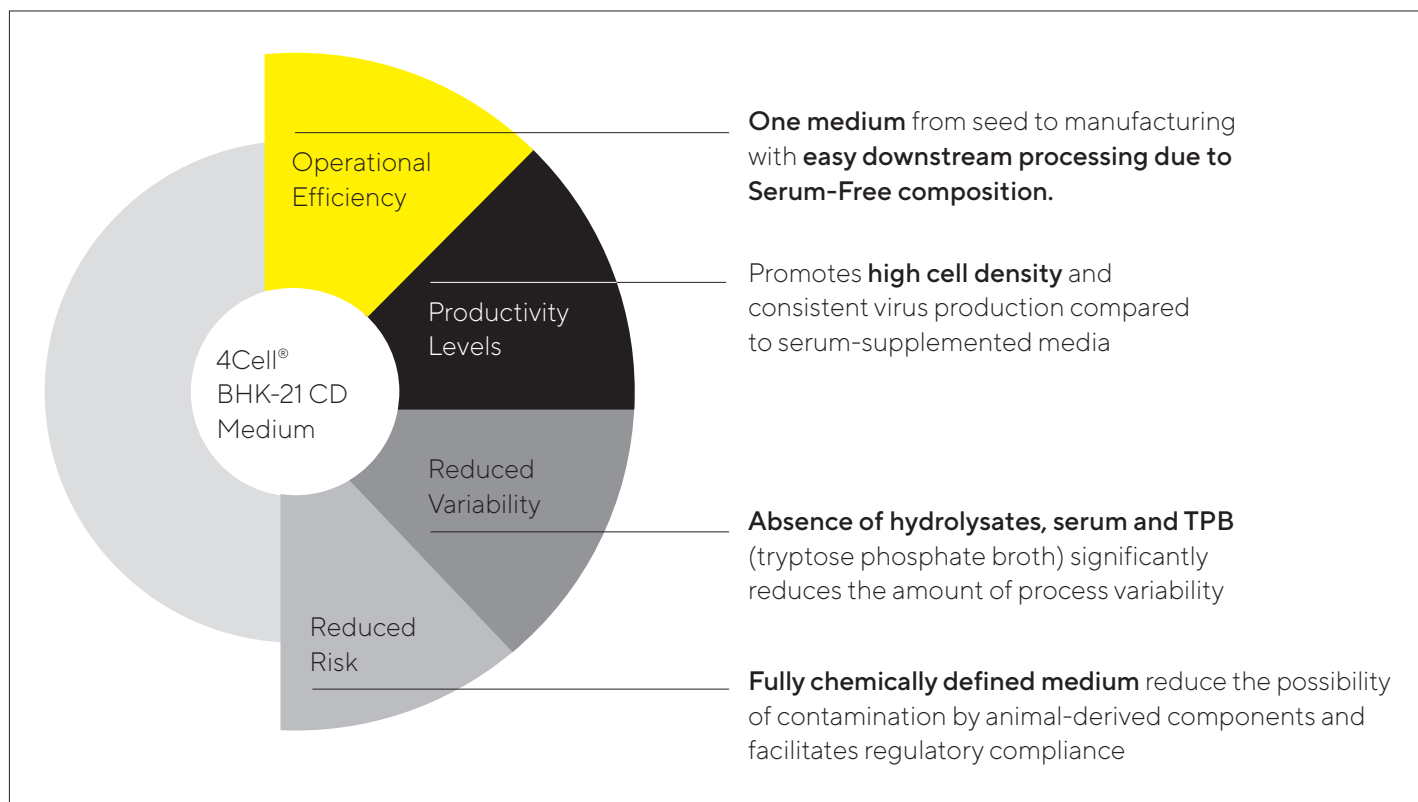
Specifications

Media Type	Chemically-defined medium: does not contain serum or undefined components
Easy-to-use	Available in liquid or complete powder
Cell Line	Developed for suspension BHK-21 (Baby Hamster Kidney) cells
Storage Condition	2-8°C, protect from light

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 further manufacturing use
- Product available in powder format
- Product available in liquid format
- AF** **Antibiotic-free:** The formulation does not contain any antibiotics

Your Benefits at a Glance



Comparison Cell Growth Performance Between BHK-21 & Reference Media

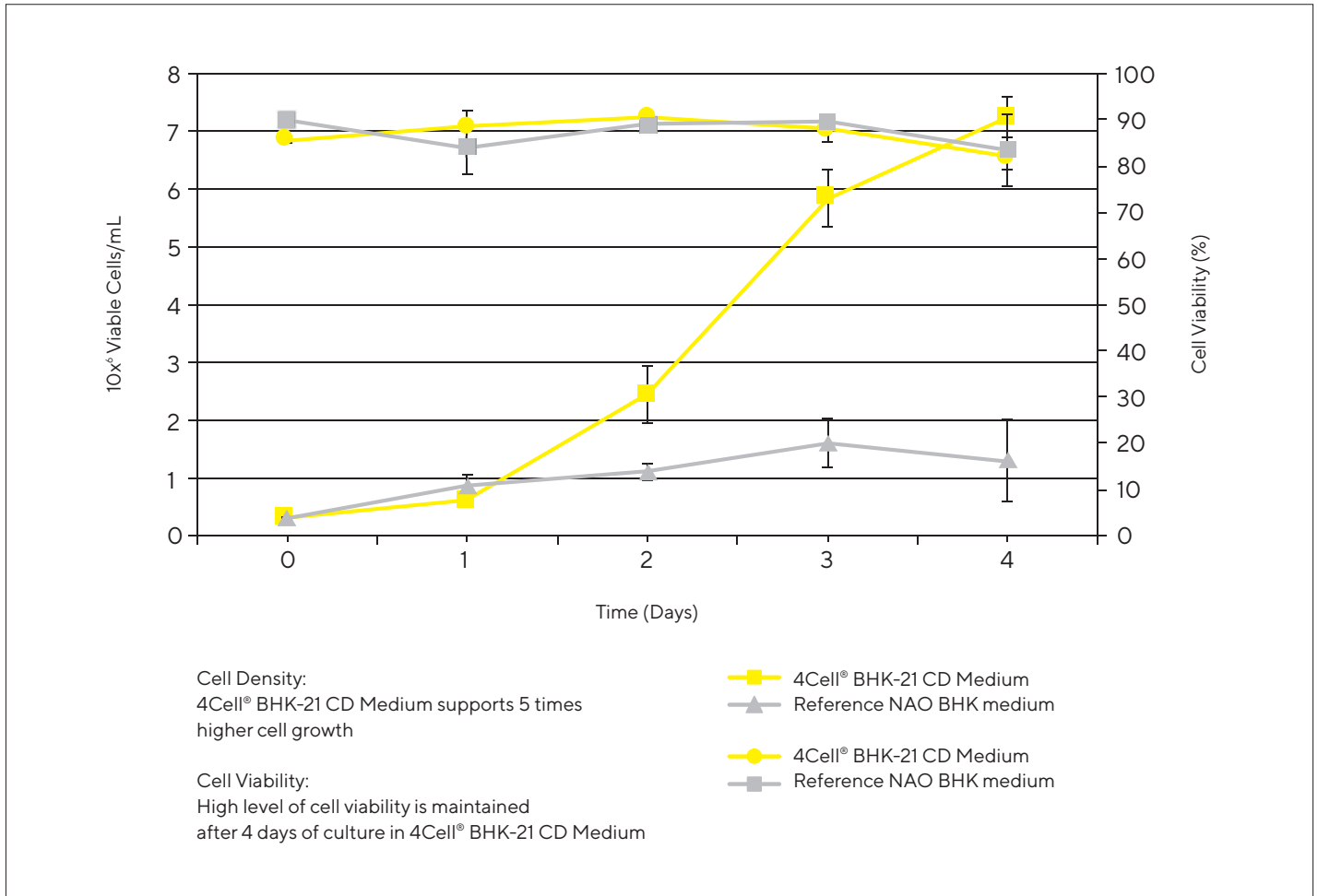


Figure 1: BHK-21 cell growth and viability in BHK-21 CD, serum-free, non-animal origin media compared to reference BHK-21 serum-free, non-animal origin medium. BHK-21 cells were sequentially adapted to 4Cell® BHK-21 CD Medium, as well as to serum-free, non-animal origin medium. Once the cells were adapted to each test medium, growth curves were determined by seeding 3×10^5 viable cells/mL in 4 L (working volume) STR bioreactors. Each test condition represents the average of 2 replicates.

Ordering Information

Description	Form	Package	Volume Size*	Order Code
4Cell® BHK-21 CD Medium	Liquid	Bottle	500 mL 6 × 500 mL Bottle	CFV3FA0001
4Cell® BHK-21 CD Medium	Liquid	Bottle	500 mL 2 × 500 mL Bottle	CFV3FA0002
4Cell® BHK-21 CD Medium	Powder	Bucket	10 L 1 Bucket	CQV3FA0010
4Cell® BHK-21 CD Medium	Powder	Bucket	50 L 1 Bucket	CQV3FA0011
4Cell® BHK-21 CD Medium	Powder	Bucket	100 L 1 Bucket	CQV3FA0014

* Other sizes are available on request

For more information please contact:
Europe | Asia: CellCultureMedia.EU@sartorius.com
Americas: CellCultureMedia.NA@sartorius.com


Or your Sartorius sales representative.

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen
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