

CHO CD EfficientFeed™ B AGT™ Nutrient Supplement

Description

CHO CD EfficientFeed™ B AGT™ Nutrient Supplement is a dry format single part supplement for fed batch addition to the culture of multiple cell lines. CHO CD EfficientFeed™ B AGT™ Nutrient Supplement is an animal origin-free, chemically defined formulation that contains no proteins, hydrolysates, or components of unknown composition.

Product	Catalog No.	Amount	Storage	Shelf Life*
CHO CD EfficientFeed™ B AGT™ Nutrient Supplement**	A12456-05 A12456-01	1 × 10 L 1 × 100 L	2°C to 8°C; Store Dark and Dry.	24 months

* Shelf Life duration is determined from Date of Manufacture.

** AGT = Advanced Granulation Technology.

Product Use

For Research or Further Manufacturing Use. Not for use in diagnostic procedures.

Safety Information

Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Prepare Medium

Reconstitute CHO CD EfficientFeed™ B AGT™ Nutrient Supplement

1. Measure room temperature (15°C to 30°C) deionized or distilled water to 90% of final volume into an appropriately sized clean vessel.
2. Add CHO CD EfficientFeed™ B AGT™ at 54.15 g/L. Mix for a minimum of 30 minutes. Solution will be cloudy, improving with time.
3. Slowly adjust to pH 7.20 with 5N Hydrochloric Acid. Mix for 15–30 minutes.
4. Add room temperature deionized or distilled water to final volume. Mix for 15 minutes.
5. Sterilize by membrane filtration.

Note: Once product is filter sterilized, use immediately or dispense product into single-use aliquots prior to storage. Store product at 2°C to 8°C (protected from light) up to 6 months.

Use

CHO CD EfficientFeed™ B AGT™ Nutrient Supplement can be added as multiple discrete additions every 1–3 days or as a continuous feed. Total feed volume can range from 10–40% of the original culture volume depending on the cell line and process.

Related Products

Product	Catalog No.
Water, Distilled	15230
Water for Injection (WFI) for Cell Culture,	A12873
CD FortiCHO™ Medium	A11483
CD OptiCHO™ Medium	12681
CD CHO Medium	10743
CHO CD EfficientFeed™ A	A10234
CHO CD EfficientFeed™ B Liquid Nutrient Supplement	A10240
CD EfficientFeed™ C AGT™	A13275
Trypan Blue Stain	15250
Countess® Automated Cell Counter	C10227

Explanation of Symbols and Warnings

The symbols present on the product label are explained below:

			
Temperature Limitation	Manufacturer	Batch code	Use By:
			
Caution, consult accompanying documents	Consult instructions for use	Sterilized using aseptic processing techniques	
			
Read Safety Data Sheet	Catalog number	Protect from light	

Limited Use Label License: Internal Research and Bioproduction Use

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product (a) to perform internal research for the sole benefit of the purchaser; and (b) to culture cells for the purpose of producing a product wherein the product will be used for any or all of the following: (i) internal research use by the purchaser; (ii) resale for internal research use by third parties; (iii) performance of research conducted by the purchaser on a fee for service or contract basis for or on behalf of third parties; (iv) resale for use as a human therapeutic agent or diagnostics product or component by third parties; (v) performance of manufacturing services conducted by the purchaser on a fee for service or contract basis for or on behalf of third parties.

The purchase of this product does not grant the purchaser any additional rights, including (without limitation) the right to transfer or resell the product in any form, the right to use the product as a therapeutic agent or diagnostics test component, or to use the product to perform other tests on a contract or fee per test basis for or on behalf of third parties. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

For additional technical information such as Safety Data Sheets (SDS), Certificates of Analysis, visit www.lifetechnologies.com/support
For further assistance, email techsupport@lifetech.com

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

www.lifetechnologies.com

