

# CHO-S™ Cells (cGMP Banked) and Media Kit

Catalog Number A11557-01

Pub. No. MAN0007378 Rev. 3.0



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

#### **Product description**

Gibco™ CHO-S™ Cells (cGMP Banked) and Media Kit have been developed for the growth of Chinese Hamster Ovary (CHO) cells and expression of recombinant proteins in suspension culture. CHO-S™ cells have been adapted to CD CHO Medium for serum–free suspension growth, and subsequently banked and tested to meet cGMP quality standards. CD CHO Medium is an animal origin-free (AOF), chemically defined medium that contains no proteins, hydrolysates, or components of unknown composition.CD CHO Medium is formulated without L-glutamine for greater stability, and without phenol red to minimize potential for estrogen-like effects. CD CHO Medium is made without hypoxanthine and thymidine for use in dihydrofolate reductase (DHFR) amplified systems.

#### Contents and storage

Table 1 CHO-S™ Cells (cGMP Banked) and Media Kit, Cat. No. A11557-01

Contents	Cat. No.	Amount	Storage	Shelf life <sup>[1]</sup>
CD CH0 Medium	10743-029	1000 mL	2°C to 8°C. Protect from light.	18 months
CHO-S™ Cells (cGMP Banked)	A11364-01	1 vial <sup>[2]</sup>	−200°C to −125°C. Liquid Nitrogen.	_
L-Glutamine, 200 mM	A2916801	100 mL	−20°C to −5°C. Protect from light.	24 months

<sup>[1]</sup> Shelf-Life duration is determined from Date of Manufacture.

#### Important information

- CHO-S<sup>™</sup> cells have been produced, banked, and tested to meet current Good Manufacturing Practice regulations 21 CFR Parts 210, 211, 600, and 610.
- CHO-S<sup>™</sup> Cells: Stable when maintained at -200°C to -125°C.

#### Prepare medium

CD CHO Medium requires supplementation with L-glutamine prior to use.

- 1. As eptically add L-glutamine to 8 mM final concentration (40 m L/L), to the medium before use.
- 2. If cell clumping occurs, add 1 mL/L of Anti-Clumping Agent to medium. After any thaw or changes in media composition, subculture cells for a minimum of 3 passages before use in other applications.

**Note:** Consider reducing L-glutamine concentration for fed batch or perfusion protocols, or to reduce ammonia levels.

**Note:** Addition of a surfactant (e.g., Pluronic  $^{\text{TM}}$  F-68) is not required.

#### Culture conditions

**Medium**: Complete CD CHO Medium **Cell line**: CHO-S<sup>™</sup> Cells (cGMP Banked)

Culture type: Suspension

Culture vessels: Shake flask or spinner bottle

**Temperature range**: 36°C to 38°C

**Incubator atmosphere**: Humidified atmosphere of 5–10% CO<sub>2</sub> in air. Ensure proper gas exchange and minimize exposure of cultures to light.

#### Recovery

- 1. Rapidly thaw (<2 minutes) frozen vial of cells in a 37°C water bath.
- Transfer the entire contents of the cryovial into a 125-mL shake flask containing 29 mL of prewarmed complete CD CHO Medium.

If thawed properly, cell density should be  $\ge 3 \times 10^5$  viable cells/mL, and viability should be  $\ge 90\%$ .



<sup>[2] 1</sup> vial contains  $\ge 1 \times 10^7$  cells/vial

- 3. Incubate at 37°C in a humidified atmosphere of 5–10% CO<sub>2</sub> in air on an orbital shaker platform rotating at 125–135 rpm. Loosen flask caps (or use vented caps) to allow for gas exchange.
- 4. Subculture cells, 2–3 days post-thaw, when viable cell density reaches  $1 \times 10^6$  cells/mL in mid-logarithmic phase of growth. Seed cultures at a density of  $3 \times 10^5$  viable cells/mL. Subculture cells a minimum of 3 passages before use in other applications.

**Note:** Do not centrifuge CHO-S<sup>™</sup> cells as they are extremely fragile upon recovery from cryopreservation.

### Scaling up CHO-S™ Cells in CD CHO Medium

CHO-S<sup>™</sup> cultures can be scaled up in spinner bottles or stirred tank bioreactors using the following guidelines.

- Determine the optimum spinner or impeller speed for your bioreactor depending on culture requirements.
- Seeding density: We recommend an optimized seeding density of 1-2 × 10<sup>5</sup> viable cells/mL.

**Note:** If the split ratio of cells to fresh media is <1:2, we recommend to spin down the cell suspension at  $100 \times g$  for 5–10 minutes, and resuspending the cell pellet in fresh complete CD CHO Medium prior to inoculating the spinner or bioreactor culture.

#### Cryopreservation

Prepare the desired quantity of cells in a tissue culture flask, harvesting in mid-log phase of growth when viable cell density reaches  $>1 \times 10^6$  cells/mL with viability >90%.

- Determine the viable cell density and calculate the required volume of cryopreservation medium to give a final viable cell density of ≥1 × 10<sup>7</sup> cells/mL.
- 2. Prepare the required volume of cryopreservation medium (90% fresh complete, and 10% DMSO) and store at 4°C until use.

**IMPORTANT!** Prepare cryopreservation medium on the day of use.

- 3. Harvest cells by centrifugation at 100 × g for 5–10 minutes. Resuspend the pellet in the pre-determined volume of 4°C cryopreservation medium.
- Dispense aliquots of this suspension into cryovials according to the manufacturer's specifications (i.e., 1.5 mL in a 2-mL cryovial).
- **5.** Achieve cryopreservation in an automated or manual controlled rate freezing apparatus following standard procedures (1°C decrease per minute).
- Transfer frozen cells to liquid nitrogen (vapor phase); storage at −200°C to −125°C is recommended.

Check viability of cryopreserved cells 24 hours after storage of vials in liquid nitrogen. See "Recovery".

#### Related products

Product	Catalog No.	
Anti-Clumping Agent	0010057AE	
FreeStyle™ MAX Reagent	16447	
FreeStyle™ MAXCHO Expression System	K9000-20	
EfficientFeed™ A+ AGT™ Supplement	A25023	
EfficientFeed™ B+ AGT™ Supplement	A25030	
EfficientFeed™C+ AGT™ Supplement	A25031	
CD CH0 AGT™ Medium	12490	
CD CHO Medium (1X), Liquid	10743	
Water, Distilled	15230	
Freedom™ CH0-S™ Kit	A13696	
Countess™ Automated Cell Counter	C10227	
Trypan Blue Stain	15250	

## **Explanation of symbols**

		LOT		REF
Temperature Limitation	Manufacturer	Batch code	Use By	Catalog Number
<u> </u>	<u>i</u>		STERILE A	
Caution, consult accompanying documents	Consult instructions for use	Keep away from light	Sterilized using aseptic processing techniques	

#### 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.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Limited Use Label License No. 335:Technology with CRO Rights: This product and its use is the subject of one or more issued and/or pending U.S. and foreign patent applications owned or controlled by Life Technologies Corporation. The purchase of this product conveys to the buyer (whether the buyer is an academic or for-profit entity) the right under the aforementioned patents and patent applications to use the purchased amount of the product and components of the product in research. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes except that buyer may transfer this product and its components to service providers for use by such service providers in research conducted for consideration solely for the benefit of the buyer, provided that each such service provider agree in writing (a) not to transfer this product, its components, or materials or information made using this product or its components to any third party, (b) to use this product, its components, and any materials or information made using this product or its components solely for research and solely for the benefit of the buyer, and (c) to return to buyer or destroy any and all remaining product, components, or materials made using this product or its components upon completion or termination of its services for buyer. The buyer may also transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California

Limited Use Label License No. 336: Transformed Host Cell Disclaimer: This product is not provided with a license to U.S. Patent No. 6,331,415, or any patents or patent applications related thereto. Each purchaser of this product should consider whether additional rights are required for the purchaser's particular application under the above patents.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

