

CTS™ GlutaMAX™ Supplement

Catalog Numbers A1286001, A4737001, A4737003

Pub. No. MAN0007389 Rev. C



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](https://www.thermofisher.com/support).

Product description

Gibco™ CTS™ GlutaMAX™ Supplement, L-alanyl-L-glutamine, is a dipeptide substitute for L-glutamine and can be used as a direct substitute for L-glutamine at equimolar concentrations in mammalian and stem cell culture with minimal or no adaptation. CTS™ GlutaMAX™ Supplement is highly soluble, heat stable, and improves growth efficiency and performance of mammalian cell culture systems. The supplement eliminates problems associated with the spontaneous breakdown of L-glutamine into ammonia during incubation. The Gibco™ CTS™ product line enables you to reduce your burden in qualifying reagents during your transition from research applications to clinical applications.

Contents and storage

Product	Catalog No.	Amount	Storage	Shelf Life ^[1]
CTS™ GlutaMAX™ Supplement (bag)	A4737001	100 mL	15°C to 30°C	36 months
	A4737003	1 L		
CTS™ GlutaMAX™ Supplement (bottle)	A1286001	100 mL	15°C to 30°C	24 months

^[1] Shelf Life duration is determined from Date of Manufacture.

Important information

- Formulated as a 200-mM L-alanyl-L-glutamine dipeptide in 0.85% NaCl.

Related products

Product	Catalog No.
L-glutamine-200mM (100X), Liquid	25030

Use


Direct equimolar substitution for L-glutamine in mammalian cell culture.

Supplement with CTS™ GlutaMAX™ Supplement

- Aseptically add the required molar concentration to culture medium.
- If a lag in growth is observed when transitioning culture from L-glutamine to CTS™ GlutaMAX™ Supplement, it is recommended to supplement the culture media with both CTS™ GlutaMAX™ Supplement and L-glutamine at a 3:1 concentration ratio. For instance, if an 8 mM final concentration is required, you should use 6 mM CTS™ GlutaMAX™ Supplement and 2 mM L-glutamine. This supplementation strategy should be applied for the first and second passages during the adaptation to the new culture conditions.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 3175 Staley Road | Grand Island, New York 14072 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0007389 C

Revision	Date	Description
C	25 November 2024	Added Cat. No. A4737003. Updated the shelf life of CTS™ GlutaMAX™ Supplement (bag) to 36 months. Updated product name to CTS™ GlutaMAX™ Supplement. The version numbering was changed from a numerical format to a letter-based format in accordance with internal document control procedures.
2.0	5 March 2021	Rebranded. Added SKU A4737001.
1.0	31 June 2012	New document.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Limited Use Label License No. 517: Internal Research and Bioproduction Use: Notice to Purchaser: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product to (a) perform internal research for the sole benefit of the purchaser; (b) manufacture protein (or other biological material) for resale; and (c) perform research or manufacturing services conducted by the purchaser on a fee for service or contract basis for or on behalf of third parties. However, the purchaser may transfer this product, its components, or materials made using this product to a third party (including contract research/manufacturing organizations), provided that each such third party agrees in writing to use such product, components, or materials solely on behalf of the purchaser, and such third party is restricted from further transferring any such product, components, or materials to any individual or entity other than the purchaser. No additional rights are granted. By purchasing this product, the purchaser agrees not to: (1) resell the product in any form; (2) use the product as a therapeutic agent or diagnostics test component; (3) reverse engineer the product or cause the product to be reverse engineered; or (4) use the product for purposes other than what is indicated in this Limited Use Label License. Life Technologies is not aware of Intellectual Property ("IP") that would be infringed by the manufacture or sale of its products. Customers are urged to perform an IP search and analysis specific to their manufacturing processes and target biologic products. Should that analysis require information about Life Technologies products, Life Technologies will provide to the customer's IP counsel the information necessary to analyze any relevant IP in a manner that protects Life Technologies proprietary information. The purchaser is responsible for obtaining all regulatory approvals necessary for any therapeutic or diagnostic use of the protein (or biological material) manufactured using this product. For information on obtaining additional rights, please contact outlicensing@thermofisher.com Licensing and Commercial Supply, Thermo Fisher Scientific, 5823 Newton Drive, Carlsbad, California 92008 USA

©2012-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.