

CTS™ GlutaMAX™-I

Catalog Numbers A4737001 and A1286001

Pub. No. MAN0007389 Rev. 2.0



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

CTS™ GlutaMAX™-I, 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™-I is highly soluble, heat stable, and improves growth efficiency and performance of mammalian cell culture systems. CTS™ GlutaMAX™-I 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™-I (bag)	A4737001	100 mL	15°C to 30°C	12 months
CTS™ GlutaMAX™-I (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.

Use

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

Supplement with CTS™ GlutaMAX™-I

- 1. Aseptically add the required molar concentration to culture medium.
- 2. If a lag in growth is observed upon conversion to CTS[™] GlutaMAX -I, we recommend that you supplement the culture media with CTS[™] GlutaMAX[™]-I and L-glutamine at a 3:1 concentration (e.g., if an 8 mM final concentration is required, use 6 mM CTS[™] GlutaMAX[™]-I and 2 mM L-glutamine) for the first and second passages during adaptation to the new culture conditions.

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 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.

Related products

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





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.

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.

Revision history: Pub. No. MAN0007389

Revision	Date	Description
2.0	05 March 2021	Rebranded. Added Sku A4737001.
1.0	2012	New document.

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 responsible for obtaining all regulatory approvals necessary for any therapeutic or diagnostic use of the protein (or biological material) ma

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

