GlutaMAX[™]-1, 200 mM (100X)

Pub. No. MAN0007318 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

GlutaMAX[™]-1, L-alanyl-L-glutamine, is a dipeptide substitute for L-glutamine. GlutaMAX[™]-1 can be used as a direct substitute for L-glutamine at equimolar concentrations in both adherent and suspension mammalian cell cultures with minimal or no adaptation. GlutaMAX[™]-1 is highly soluble, heat-stable, and improves growth efficiency and performance of mammalian cell culture systems. GlutaMAX[™]-1 eliminates problems associated with the spontaneous breakdown of L-glutamine into ammonia during incubation, allowing for longer lasting cultures. GlutaMAX[™]-1 supplement is also available pre-supplemented in a variety of media formulations.

Product	Cat. No.	Amount	Storage	Shelf Life ^[1]
GlutaMAX™-1, 200 mM (100X)	35050061	100 mL	15°C to 30°C	24 months
	35050079	20 x 100 mL		
	35050038 ^[2]	100 mL		
	35050087 ^[2]	20 x 100 mL		

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

Important information

 Formulated as a 200 mM (100X) L-alanyl-L-glutamine dipeptide in 0.85% NaCl

Use

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

Supplement with GlutaMAX™-1

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

Related products

Product	Cat. No.			
L-glutamine 200mM (100X), liquid	25030			
GlutaMAX™-1 CTS™	A12860			
DMEM/F-12, GlutaMAX™	10565			
RPMI, GlutaMAX™	61870			
OptiMEM™ Reduced Serum Medium, GlutaMAX™	51985			

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.

thermofisher.com/support | thermofisher.com/askaquestion



Life Technologies Corporation | 3175 Staley Road | Grand Island, NY 14072 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.

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

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. For Research Use or Further Manufacturing. Not for diagnostic use or direct administration

For Research Use or Further Manufacturing. Not for diagnostic use or direct administration into humans or animals.



^[2] For European customers only.