

Recombinant Human Hepatocyte Growth Factor (HGF)

Catalog Numbers PHG0324 (10 µg), PHG0321 (100 µg), PHG0254 (10 µg),

Pub. No. MAN0004448 Rev. B.0








Product specifications

Lot number	See product label.
Molecular weight	70 kDa, single chain, glycosylated
Purity	>95% as determined by SDS PAGE analysis.
Biological activity	ED ₅₀ ≤20 ng/mL, determined by the dose-dependent stimulation of the proliferation of the monkey epithelial cell line 4MBR-5. Determine the optimal concentration for each specific application using an initial dose response assay.
Formulation	Lyophilized, with PBS.
Sterility	Filtered before lyophilization through a 0.2 micron sterile filter.
Endotoxin	<0.15 EU/µg
Production	Produced in human cell expression system and purified via sequential chromatography.
Reconstitution recommendation	Centrifuge the vial briefly, before opening to bring the contents to the bottom. Reconstitute the lyophilized protein in sterile PBS containing 0.1% human serum albumin. Apportion the reconstituted protein into working aliquots and store at ≤ -20°C. Make any further dilutions of the reconstituted protein in low endotoxin medium or a buffered solution containing a carrier protein such as heat inactivated FCS or tissue culture grade BSA.
Suggested working dilutions	The optimal concentration should be determined for each specific application.
Storage	Store the lyophilized protein at ≤ -20°C, preferably desiccated. Upon reconstitution, apportion into working aliquots and store at ≤ -20°C. Avoid repeated freeze-thaw cycles.
Expiration date	Expires one year from date of receipt when stored as instructed.
References	Rubin JS, Chan AM, Bottaro DP, Burgess WH, Taylor WG, Cech AC, Hirschfield DW, Wong J, Miki T, Finch PW, et al. (1988) A broad-spectrum human lung fibroblast-derived mitogen is a variant of hepatocyte growth factor. Proc Natl Acad Sci U S A. 1991 Jan 15;88(2):415-9. Jiang WG, Martin TA, Parr C, Davies G, Matsumoto K, Nakamura T. (2005) Hepatocyte growth factor, its receptor, and their potential value in cancer therapies. Crit Rev Oncol Hematol. 2005 Jan;53(1):35-69.

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.

Explanation of Symbols

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

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: 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.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.