

Recombinant Human Growth Hormone (HGH)

Catalog Numbers PHH0014

Pub. No. MAN0004449 Rev. 2.0








Product specifications

Amount	10 µg
Lot number	See product label.
Molecular weight	22 kDa, monomer, non-glycosylated.
Purity	≥95% as determined by SDS PAGE analysis.
Biological activity	ED ₅₀ = 0.4–2 ng/mL, determined by the dose-dependent stimulation of the proliferation of rat lymphoma line Nb2-11 cells (prolactin indicator cell line).
Formulation	Lyophilized in PBS.
Endotoxin	<1 EU/µg
Production	Recombinant human growth hormone (HGH) is expressed in human cell expression system and purified via sequential chromatography. Manufactured using all non-animal reagents, and containing only components of human or humanized origin.
Reconstitution recommendation	Centrifuge the vial briefly, prior to opening to bring the contents to the bottom. Reconstitute lyophilized protein in sterile PBS containing 0.1% human serum albumin. Make any further dilutions 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	Gout PW, Beer CT, Noble RL. (1980) Prolactin-stimulated growth of cell cultures established from malignant Nb rat lymphomas. Cancer Res. 1980 Jul;40(7):2433-6. Goffin V, Shiverick KT, Kelly PA, Martial JA. (1996) Sequence-function relationships within the expanding family of prolactin, growth hormone, placental lactogen, and related proteins in mammals. Endocr Rev. 1996 Aug;17(4):385-410.

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.

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		

Manufacturer's address: Life Technologies Corporation | 7335 Executive Way | Frederick, MD 21704 | USA

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.

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.

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

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.