

Recombinant Human Epidermal Growth Factor (EGF)

Catalog Numbers PHG0314 (10 µg), PHG0315 (25 µg), PHG0311 (100 µg), PHG0313 (1 mg)

Pub. No. MAN0003587 Rev. B.0








Product specifications

Lot number	See product label.
Molecular weight	6.2 kDa
Purity	>95% as determined by SDS PAGE analysis.
Biological activity	ED ₅₀ <0.400 ng/mL, determined by the dose dependent proliferation of mouse BALB/3T3 cells. Determine the optimal concentration for each specific application using an initial dose response assay.
Formulation	Lyophilized, carrier free.
Sterility	The protein is eluted in acetonitrile and then lyophilized under aseptic conditions.
Endotoxin	<0.1 ng/µg
Production	Produced in <i>E. coli</i> 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 buffer to a concentration of 0.1–1.0 mg/mL. 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 2°C to 8°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	<p>Gregory, H. (1975) Isolation and structure of urogastrone and its relationship to epidermal growth factor. <i>Nature</i> 257(5524):325–327.</p> <p>Ediger, TL, and Toews, ML. (2000) Synergistic stimulation of airway smooth muscle cell mitogenesis. <i>J. Pharmacol. Exp. Therapeut.</i> 294:1076–1082.</p> <p>French, JD, Tschumper RC, and Jelinek DF. (2002) Analysis of IL-6-mediated growth control of myeloma cells using a gp130 chimeric receptor approach. <i>Leukemia</i> 16(6):1189–1196.</p> <p>French, JD, Walters, DK, and Jelinek DF. (2003) Transactivation of gp130 in myeloma cells. <i>J. Immunol.</i> 170(7):3717–3723.</p> <p>Liesveld, JL, Rosell, K., Panoskaltis, N., Belanger, T., Harbol, A., and Abboud, CN. (2001) Response of human CD34(+) cells to CXC, CC, and CX3C chemokines: Implications for cell migration and activation. <i>J. Hematotherapy and Stem Cell Research</i> 10(5):643–655.</p> <p>Stephens, JM, Lumpkin, SJ, and Fishman, JB. (1998) Activation of signal transducers and activators of transcription 1 and 3 by leukemia inhibitory factor, oncostatin-M, and interferon-gamma in adipocytes. <i>J. Biol. Chem.</i> 273(47):31,408–31,416.</p>

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		

NOTE TO WRITER: ADD CONKEYREF SOURCE TO PUBLICATION

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.

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

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