

Recombinant Human Keratinocyte Growth Factor-2

Publication Number MAN0004470

Rev. 1.00

Catalog Number:	PHC1536
Quantity:	25 µg
Lot Number:	See product label.
Molecular Weight:	19.3 kDa (170 amino acid residues)
Purity:	> 95% as determined by SDS-PAGE and RP-HPLC.
Biological Activity:	ED ₅₀ <0.5 ng/mL (corresponding to a specific activity of >2 × 10 ⁶ units/mg). The biological activity was determined by measuring the dose dependent incorporation of ³ H-thymidine by BaF3 cells. The optimal concentration for each specific application should be determined by an initial dose-response assay.
Formulation:	Lyophilized from a solution containing 1 mg/mL recombinant Keratinocyte Growth Factor-2, 5 mM sodium phosphate, pH 7.4, and 80 mM NaCl.
Sterility:	Sterile-filtered prior to lyophilization.
Endotoxin:	<0.1 ng/µg
Production:	Recombinant human Keratinocyte Growth Factor-2 is produced in <i>E. coli</i> and purified via sequential chromatography.
Reconstitution Recommendation:	We recommend that the vial be briefly centrifuged prior to opening to bring the contents to the bottom. Reconstitute lyophilized recombinant human Keratinocyte Growth Factor-2 in sterile, distilled water to a concentration greater than 0.1 mg/mL. Further dilutions should be made in buffered solution containing a carrier protein, such as 0.1% HSA or BSA.
Suggested Working Dilutions:	The optimal concentration should be determined for each specific application.
Storage:	This lyophilized protein is stable at room temperature for three weeks, but should be kept at -20°C for long term storage. Upon reconstitution, KGF-2 can be stored at 4°C for 2-7 days. For long term storage, apportion recombinant human KGF-2 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:	Steiling, H. and S. Werner (2003) Fibroblast growth factors: key players in epithelial morphogenesis, repair and cytoprotection. <i>Curr. Opin. Biotechnol.</i> 14:533-537. Hogan, B.L. (1999) Morphogenesis. <i>Cell</i> 96:225-233. Werner, S. (1998) Keratinocyte growth factor: a unique player in epithelial repair processes. <i>Cytokine Growth Factor Rev.</i> 9:153-165.

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description
	Catalog Number
	Research Use Only
	Use by
	Manufacturer
	Without, does not contain
	Protect from light
	Directs the user to consult instructions for use (IFU), accompanying the product.

Symbol	Description
	Batch code
	In vitro diagnostic medical device
	Temperature limitation
	European Community authorized representative
	With, contains
	Consult accompanying documents

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

For Research Use Only. Caution: Not for human or animal therapeutic or diagnostic use.

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

Revision Date 01 June 2011

