

Recombinant Rat GRO








Publication Number MAN0004497







Rev. 1.00

Catalog Number:	PRC1066
Quantity:	25 µg
Lot Number:	See product label.
Molecular Weight:	Approximately 7.8 kDa, 72 amino acid residues.
Purity:	Recombinant rat GRO is >98% pure by SDS-PAGE and HPLC analyses.
Biological Activity:	The biological activity of this recombinant rat GRO is determined by measuring concentration dependent chemotaxis with isolated rat neutrophils. Significant chemotaxis is observed at a concentration range of 10-100 ng/mL. The optimal concentration should be determined for each specific application. This protein is also known as GRO alpha, KC, and CINC.
Formulation:	Lyophilized, carrier-free, preservative-free. Sterile filtered prior to lyophilization.
Endotoxin:	<0.1 ng/µg (1 EU/µg) of rat GRO.
Production:	Recombinant rat GRO is produced in <i>E. coli</i> .
Reconstitution Recommendation:	We recommend that this vial be briefly centrifuged prior to opening to bring the contents to the bottom. Recombinant rat GRO should be reconstituted in sterile, distilled water to a concentration of 0.1–1.0 mg/mL. Recombinant rat GRO is also soluble in most buffered solutions. Further dilutions should be made in buffered solution with carrier protein such as PBS with 1% BSA.
Storage:	This lyophilized preparation is stable at 2°C to 8°C, but should be kept at –20°C for long term storage, preferably desiccated. Upon reconstitution, the preparation is stable for up to one week at 2°C to 8°C. For maximal stability, apportion the reconstituted preparation 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:	Fekete, Z., C.J. Hauser, J.M. Adams, C.A. Adams Jr., R.M. Forsythe, G. Hasko, D.Z. Xu, D.H. Livingston, and E.A. Deitch (2001) Injury-enhanced calcium mobilization in circulating rat neutrophils models human PMN responses. Shock 16(1):15–20.

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 03 June 2011

