

Recombinant Rat Interleukin-1 (rIL-1)








Publication Number MAN0004507





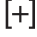

Rev. 1.00

Catalog Number:	PRC0014	PRC0015	PRC0013
Quantity:	10 µg	25 µg	1 mg
Lot Number:	See product label.		
Molecular Weight:	18 kDa		
Purity:	≥95% as determined by SDS-PAGE analysis.		
Biological Activity:	ED ₅₀ range = ≤0.005 ng/mL, determined by the dose dependent proliferation of mouse D10S cells. The optimal concentration for individual application should be determined by an initial dose response assay.		
Formulation:	Lyophilized, carrier-free.		
Sterility:	Filtered prior to lyophilization through a 0.22 micron sterile filter.		
Endotoxin:	<0.1 ng/µg		
Production:	Produced in <i>E. coli</i> and purified by sequential chromatography.		
Reconstitution Recommendation:	We recommend that this vial be briefly centrifuged prior to opening to bring the contents to the bottom. Reconstitute lyophilized rIL-1α in sterile, distilled water to a concentration of 0.1–1.0 mg/mL. Recombinant rIL-1α is also soluble in most buffered solutions. Further dilutions should be made in buffered solution containing a carrier protein such as PBS + 0.1% BSA.		
Suggested Working Dilutions:	The optimal concentration should be determined for each specific application.		
Storage:	Store lyophilized rIL-1α 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:	See product label.		
References:	<p>Nishida, T., N. Nishino, M. Takano, Y. Sekiguchi, K. Kawai, K. Mizuno, S. Nakai, Y. Masui, and Y. Hirai (1989) Molecular cloning and expression of rat interleukin-1 alpha cDNA. <i>J. Biochem.</i> 105:351–357.</p> <p>Orencole, S.F. and C.A. Dinarello (1989) Characterization of a subclone (D10S) of the D10.G4.1 helper T-cell line which proliferates to attomolar concentrations of interleukin-1 in the absence of mitogens. <i>Cytokine</i> 1:14–22.</p> <p>Hirano, T., (1994) In: <i>The Cytokine Handbook</i>, 2nd Edition, (ed A. Thomson), Academic Press Limited, San Diego, Ca, pp. 145–168.</p> <p>Dinarello, C. (1996) Biological basis for interleukin-1 in disease. <i>Blood</i> 87:2095–2147.</p> <p>Anforth, H.R., R. Bluthe, A. Bristow, S. Hopkins, M. Lenczowski, G. Luheshi, J. Lundkvist, B. Michaud, Y. Mistry, A. Van Dam, C. Zhen, R. Dantzer, S. Poole, N. Rothwell, F. Tilders, and E. Wollman (1998) Biological activity and brain actions of recombinant rat interleukin-1alpha and interleukin-1beta. <i>Eur. Cytokine Netw.</i> 9:278–288.</p>		

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.

Manufactured under ISO 13485 Quality Standard

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

