

Recombinant Mouse Macrophage Inflammatory Protein–1 Beta (MIP–1β)








Publication Part Number MAN0004605





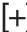

Rev. 1.00

Catalog Number:	PMC1034	PMC1031
Quantity:	10 µg	100 µg
Lot Number:	See product label.	
Molecular Weight:	~8 kD	
Purity:	≥95% pure by SDS–PAGE	
Biological Activity:	The biological activity was determined by measuring the dose dependent mobilization of intracellular calcium (calcium flux) with K562 cells transfected with human CCR5. Significant calcium mobilization is observed with ≥500 ng/mL of recombinant mouse MIP–1β. The optimal concentration for each specific application should be determined by an initial dose response assay.	
Formulation:	Lyophilized, carrier-free.	
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. Lyophilized mouse MIP–1β should be reconstituted in sterile deionized water to a concentration of 0.1–1.0 mg/mL to regain full activity. Stock solutions should be apportioned into working aliquots and stored at ≤ –20°C. Further dilutions should be made in medium or buffered solution containing carrier protein, such as PBS with 0.1% BSA.	
Suggested Working Dilutions:	The optimal concentration should be determined for each specific application.	
Storage:	Lyophilized mouse MIP–1β should be stored at 2°C to 8°C, preferably desiccated. Store reconstituted mouse MIP–1β at ≤ –20°C (not in a frost-free freezer). Keep freeze-thaw cycles to a minimum.	
Expiration Date:	Expires one year from date of receipt when stored as instructed.	
References:	<p>Sherry, B., P. Tekamp-Olson, C. Gallegos, D. Bauer, G. Davatelis, S.D. Wolpe, F. Masiarz, D. Coit, and A. Cerami (1988) Resolution of the two components of macrophage inflammatory protein 1, and cloning and characterization of one of those components, macrophage inflammatory protein 1 beta. <i>J. Exp. Med.</i> 168:2251–2259.</p> <p>Broxmeyer, H.E., B. Sherry, S. Cooper, F.W. Ruscetti, D.E. Williams, P. Arosio, B.S. Kwon, and A. Cerami (1991) Macrophage inflammatory protein (MIP)-1 beta abrogates the capacity of MIP–1 alpha to suppress myeloid progenitor cell growth. <i>J. Immunol.</i> 147:2586–2594.</p> <p>Driscoll, K.E. (1994) Macrophage inflammatory proteins: Biology and role in pulmonary inflammation. <i>Exp. Lung Res.</i> 20:473–490.</p> <p>Ganju, R.K., P. Dutt, L. Wu, W. Newman, H. Avraham, S. Avraham, and J.E. Groopman (1998) Beta-chemokine receptor CCR5 signals via the novel tyrosine kinase RAFTK. <i>Blood</i> 91:791–797.</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 12 July 2011

