

Recombinant Human Interleukin–5 (IL–5)

Publication Number MAN0004231

Revision Date 04 May 2011

Catalog Number:	PHC0055			
Quantity:	10 µg			
Lot Number:	See product label.			
Molecular Weight:	26.0 kDa, disulfide linked homodimer; each subunit contains 116 amino acid residues.			
Description:	IL–5 is produced by T cells, mast cells, and eosinophils. IL–5 stimulates the growth and differentiation of B cells and eosinophils.			
Source:	Recombinant human IL-5 is produced in <i>E. coli</i> and purified by sequential chromatography.			
Purity:	Recombinant human IL–5 is ≥98% pure as determined by SDS-PAGE and HPLC analyses.			
Biological Activity:	The expected ED_{50} is ≤ 0.15 ng/mL (specific activity $\geq 6.6 \times 10^6$ units/mg). The biological activity is determined by measuring the IL-5 dose dependent proliferation of TF-1 cells. The optimal concentration should be determined for each specific application.			
Formulation:	Lyophilized in 20 mM Sodium Phosphate, pH 7.5. Sterile filtered prior to lyophilization.			
Reconstitution Recommendation:	We recommend that the vial be briefly centrifuged prior to opening to bring the contents to the bottom. Reconstitute recombinant human IL–5 in sterile distilled water to a concentration of 0.1–1.0 mg/mL. <i>Do not vortex</i> . This solution can be stored at 2°C to 8°C for up to 1 week. For extended storage, it is recommended to further dilute in a buffer containing a carrier protein, such as 0.1% BSA and store in working aliquots at –20°C to –80°C.			
Storage:	The lyophilized protein is stable at room temperature for 1 month, but should be kept at -20°C for long term storage. Working aliquots stored with a carrier protein are stable for at least 12 months at -20°C to -80°C. Avoid repeated freeze/thaw cycles.			
Expiration Date:	See product label.			
References:	Gubina, E., X. Luo, E. Kwon, K. Sakamoto, Y.F. Shi, and R.A. Mufson (2001) βc cytokine receptor-induced stimulation of cAMP response element binding protein phosphorylation requires protein kinase C in myeloid cells: A novel cytokine signal transduction cascade. J. Immunol. 167(8):4303–4310.			

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description	Symbol	Description
REF	Catalog Number	LOT	Batch code
RUO	Research Use Only	IVD	In vitro diagnostic medical device
Σ	Use by	X	Temperature limitation
	Manufacturer	EC REP	European Community authorized representative
[-]	Without, does not contain	[+]	With, contains
erolac, from Light	Protect from light	Â	Consult accompanying documents
ĺÌ	Directs the user to consult instructions for use (IFU), accompanying the product.		

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 <u>outlicensing@lifetech.com</u> 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

