



Certificate of Analysis

Human Interleukin-3 (Multi-Colony Stimulating Factor)

Catalog No: 11815-1

Lot No: 5888

Size: 10 µg/vial

Expiration: October 11, 2014

Description: Recombinant Human Interleukin-3

Source: A DNA sequence encoding Human IL-3 protein (Yang, Y.C. *et al.*, 1986 *Cell* 47:10) was expressed in *E. coli*

Reconstitution: Sterile PBS containing at least 0.1% human serum albumin or bovine serum albumin to prepare a stock of no less than 100 µg/ml of the cytokine.

Buffer: lyophilized in PBS containing 50 µg of bovine serum albumin per 1 µg of cytokine

Molecular Weight: Approximately 15 kDa

Purity: ≥ 97%

Activity: ED₅₀ is typically 0.1-0.4 ng/ml

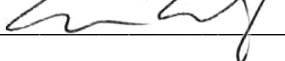
Assay Used to Measure Bioactivity: Cell proliferation assay using a human, factor -dependent cell line TF-1 (Kitamura, T. *et al.*, 1989, *J. Cell Physiol.* 140:323-334).

Shipping Conditions: Wet ice

Physical State of Product During Shipping: Lyophilized

Storage Conditions/Comments: After receipt, this product (as supplied) should be kept at -20 to -70°C for up to 1 year for retention of full activity. After reconstitution, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Under sterile conditions, this product may be kept at 2 to 8°C for up to 1 month and -20 to -70°C for up to 3 months after reconstitution. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____ 

Date: October 11, 2013

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.