



6405 Mira Mesa Blvd, Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Mouse Interleukin-3 Recombinant

Catalog #: 90189
Lot#: 240304

Description: Recombinant mouse IL-3 is a disulfide-linked monomer consisting of 135 amino acid residues and migrates as an approximately 15 kDa protein under non-reducing conditions and reducing conditions in SDS-PAGE.

Source: Optimized DNA sequence encoding murine IL-3 mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 µm filtered PBS solution.

Reconstitution: A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8°C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >98%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/µg (1 EU/µg), using the LAL gel clot method.

Biological Activity: The ED50, determined by dose-dependent stimulation of the proliferation of mouse M-NFS-60 cells, was found to be <0.05 ng/ml.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com