



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

## Human Interleukin-4 Recombinant

Catalog #: 90193-A  
Lot#: 220722

**Description:** Recombinant human interleukin-4 (IL-4), also known as B-Cell Growth Factor (BCGF), is a disulfide-linked monomeric protein consisting of 130 amino acid residue subunits and migrates as an approximately 15 kDa protein under reducing and non-reducing conditions.

**Source:** Optimized DNA sequence encoding human IL-4 mature chain was expressed in *E. coli*.

**Formulation:** Lyophilized from a 0.2 µm filtered solution containing 2.5% glycine, 0.5% sucrose, 0.01% Tween-80, 5 mM glutamic acid, pH 4.5.

**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:** >95%, 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 the dose-dependent stimulation of the proliferation of human TF-1 cells, is ≤ 0.1 ng/ml, corresponding to a specific activity of > 1x 10<sup>7</sup> units/mg.

### **Protein Sequence:**

MGLTSQLLPPLFFLLACAGNFVHGHKCD  
ITLQEIIKTLNSLTEQKTLCTELTVTDIFAA  
SKNTTEKETFCRAATVLRQFYSHHEKDT  
RCLGATAQQFHRHKQLIRFLKRLDRNL  
WGLAGLNSCPVKEANQSTLENFLERL  
KTIMREKYSK CSS

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](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)