

IL-3 (human), (recombinant)

Interleukin-3 (IL-3) is a cytokine produced by activated T cells and mast cells. IL-3 is able to induce the differentiation of hematopoietic stem cells to precursor cells of myeloid lineage (erythrocytes, megakaryocytes, granulocytes, dendritic cells and monocytes). IL-3 also has functions in the nervous system and appears to be important in several chronic inflammatory diseases.

Ordering Information

[Order Online »](#)

ALX-201-820-0010	10µg
ALX-201-820-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

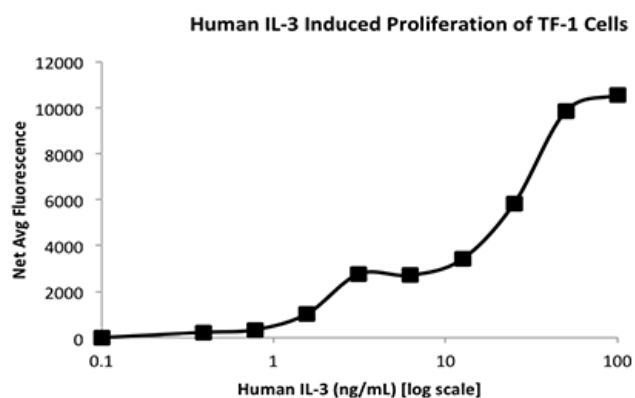


Figure 2: TF-1 cells were cultured with 0 to 100 ng/mL Human IL-3. Cell proliferation was measured after 72 hours and the linear portion of the curve can be used to calculate the ED50.

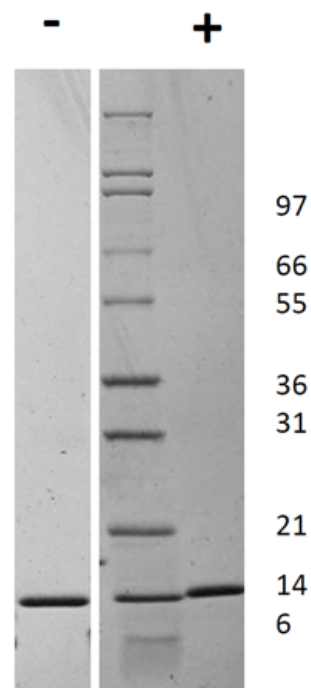


Figure 1: 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human IL-3 has a predicted MW of 15 kDa.

Handling & Storage

Use/Stability	It is recommended that a carrier protein (0.1% BSA) is added for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin-3, MCGF, Multi-CSF, HCGF, P-cell stimulation factor
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the dose-dependent stimulation of TF-1 cells which is typically observed at concentrations ED50 ≤2 ng/mL.
Endotoxin Content	≤1 EU/μg protein measured by kinetic LAL analysis.
Formulation	Lyophilized. Sterile filtered.
MW	~15kDa
Purity	≥95% (Reducing and Non-reducing SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 134 amino acids.
UniProt ID	P08700