

# 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

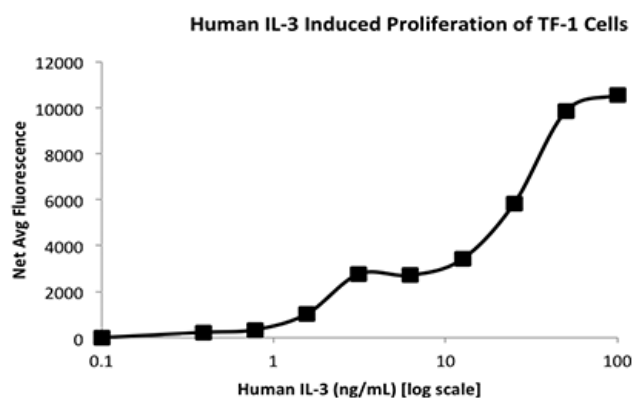
[Order Online »](#)

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

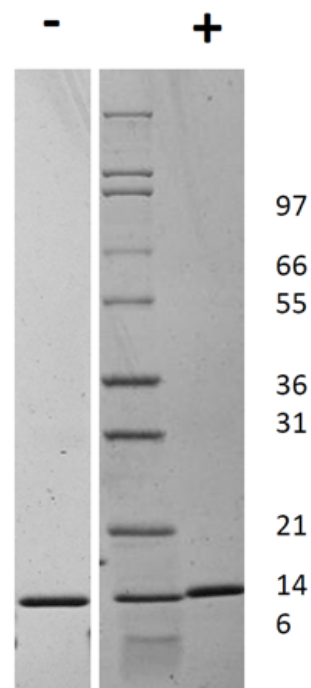
## Manuals, SDS & CofA

[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

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

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Interleukin-3, MCGF, Multi-CSF, HCGF, P-cell stimulation factor
<b>Appearance</b>	White lyophilized (freeze-dried) powder.
<b>Biological Activity</b>	The activity is determined by the dose-dependent stimulation of TF-1 cells which is typically observed at concentrations $ED_{50} \leq 2$ ng/mL.
<b>Endotoxin Content</b>	$\leq 1$ EU/ $\mu$ g protein measured by kinetic LAL analysis.
<b>Formulation</b>	Lyophilized. Sterile filtered.
<b>MW</b>	~15kDa
<b>Purity</b>	$\geq 95\%$ (Reducing and Non-reducing SDS-PAGE)
<b>Reconstitution</b>	Reconstitute in sterile water at 0.1 mg/mL.
<b>Source</b>	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 134 amino acids.
<b>UniProt ID</b>	P08700



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)