

# IL-3 (mouse), (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

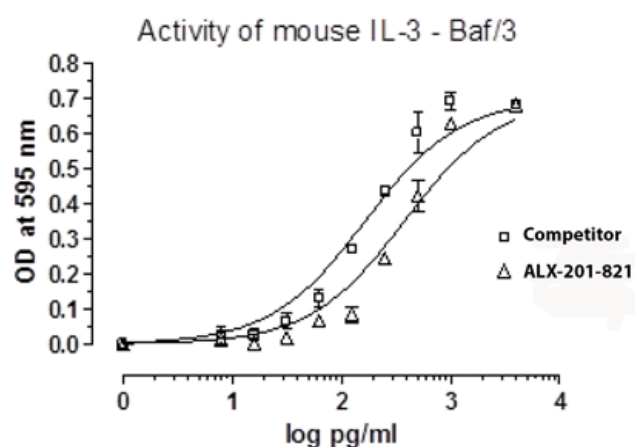
[Order Online »](#)

ALX-201-821-0010	10µg
ALX-201-821-0100	100µg

## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 1: Mouse IL-3 Bioactivity Data.** Serial dilutions of mouse IL-3, starting at 0.350ng/ml, were added to Baf/3 cells. After 41 hours cell proliferation was measured and the linear portion of the curve was used to determine the ED50. The ED50 for this lot of mouse IL-3 was 0.3-0.47ng/ml. This value is comparable to the expected range of less than 0.5ng/ml.

## Handling & Storage

<b>Use/Stability</b>	It is recommended that a carrier protein (0.1% HSA or 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 <0.1 ng/ml.
<b>Endotoxin Content</b>	≤1 EU/μg using EU/μg using USP/EP 2.6.14.
<b>Formulation</b>	Lyophilized. From sterile filtered solution containing 0.1% TFA.
<b>MW</b>	~15.1 kDa
<b>Purity</b>	≥95% (reducing and non-reducing SDS-PAGE)
<b>Reconstitution</b>	When reconstituting the product, gently pipet and wash down the sides of the vial to ensure full recovery of the protein into solution. It is recommended to reconstitute the lyophilized product with sterile water at a concentration of 0.1 mg/ml, which can be further diluted into other aqueous solutions.
<b>Source</b>	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 135 amino acids.
<b>UniProt ID</b>	P01586



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)