

# IL-6 (mouse), (recombinant)

Interleukin 6 (IL-6) is an important pro-inflammatory and anti-inflammatory cytokine expressed by T cells, macrophages and muscle cells. IL-6 signals through a receptor complex containing two receptors, IL-6R $\alpha$  and gp130. IL-6 has an important function in promoting fever and can serve to stimulate an immune response to trauma. IL-6 is often used for growth of hybridoma cell lines. Human IL-6 is active on mouse and rat cells.

## Ordering Information

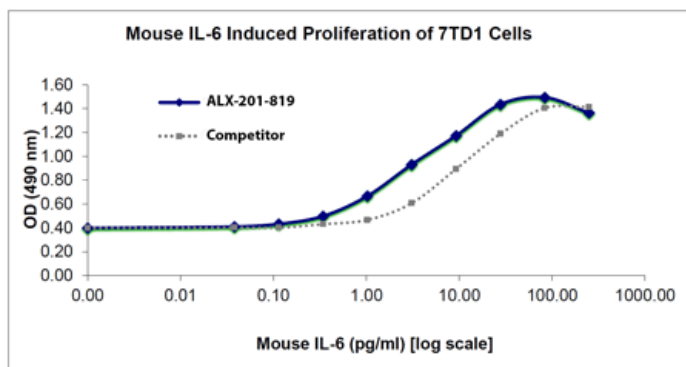
[Order Online »](#)

ALX-201-819-0010	10 $\mu$ g
ALX-201-819-0100	100 $\mu$ g

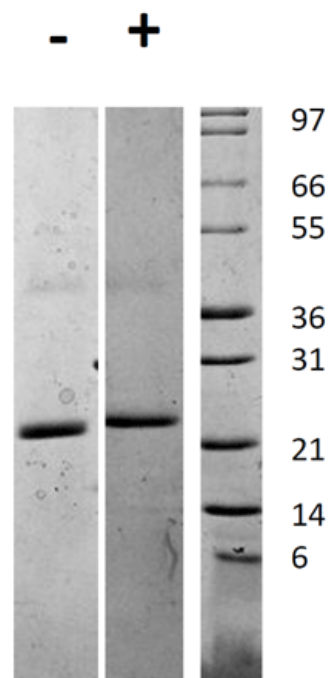
## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 1: Mouse IL-6 Bioactivity Data.** 7TD1 cells were cultured with 0 to 250pg/ml mouse IL-6. Cell proliferation was measured after 66 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of mouse IL-6 was 2.4-3.6pg/ml. This value is comparable to expected range of less than 20pg/ml.



**Figure 2:** 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Mouse IL-6 has a predicted MW of 21.7 kDa.

## 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-6, IFN- $\beta$ 2, B-cell differentiation factor, BSF-2, HSF, MGI-2
<b>Appearance</b>	White lyophilized (freeze-dried) powder.
<b>Biological Activity</b>	The activity is determined by dose-dependent induction of 7TD1 cell proliferation which is typically observed at concentrations <0.02ng/ml.
<b>Endotoxin Content</b>	≤1 EU/μg protein measured by kinetic LAL analysis.
<b>Formulation</b>	Lyophilized. Sterile filtered.
<b>MW</b>	~21.7kDa
<b>Purity</b>	≥95% (HPLC, Reducing and Non-reducing SDS-PAGE, UV spectroscopy at 280 nm)
<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 mouse IL-6 in 10 mM HCl to 0.1-1.0 mg/ml, which can be further diluted to other aqueous solutions.
<b>Source</b>	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 187 amino acids.
<b>UniProt ID</b>	P08505



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)