

# IL-6 (human), (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.

Citations: 2

[View Online »](#)

## Ordering Information

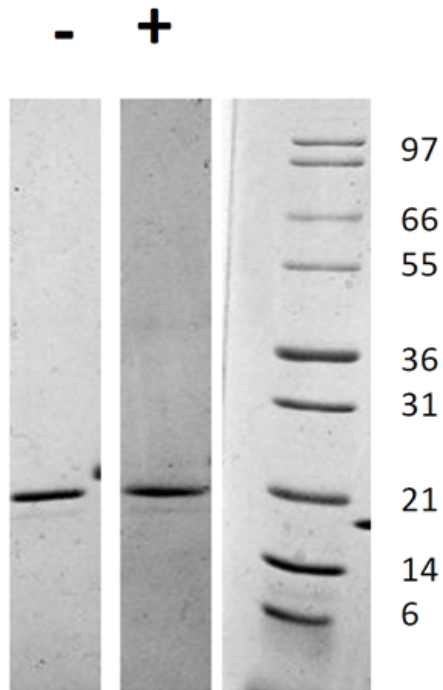
[Order Online »](#)

ALX-201-818-0020	20 $\mu$ g
ALX-201-818-0100	100 $\mu$ g

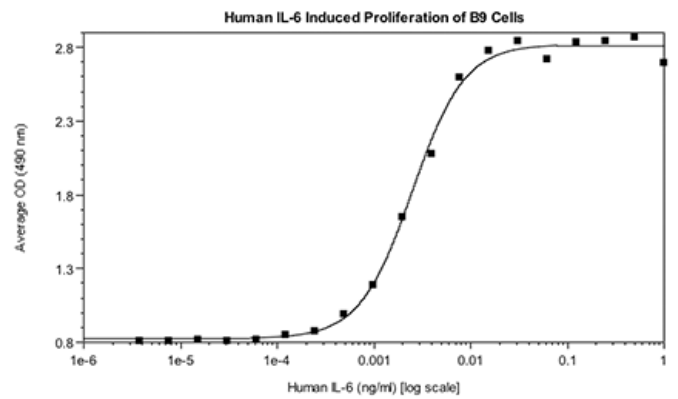
## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 1:** 1 $\mu$ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human IL-6 has a predicted MW of 20.9 kDa.



**Figure 2:** B9 cells were cultured with dilutions of Human IL-6. Cell proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

## Handling & Storage

**Use/Stability** Add 0.1% BSA for long term storage.

**Handling** Store at -80°C after reconstitution

**Long Term Storage** -80°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Interleukin-6, IFN-β2, B-cell differentiation factor, BSF-2, HSF, MGI-2

**Appearance** White lyophilized (freeze-dried) powder.

**Biological Activity** ≤0.025 ng/mL. The ED50 determined by B9 cell proliferation assay

**Endotoxin Content** ≤1 EU/μg using USP /EP 2.6.14

**Formulation** Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 10mM sodium phosphate, pH 7.5.

**MW** ~21kDa

**Purity** ≥97% using Coomassie Staining

**Reconstitution** Sterile water at 0.1 mg/mL

**Source** Produced in *E. coli*. Non-glycosylated protein, containing 184 amino acids.

**UniProt ID** P05231

**Last modified: May 29, 2024**



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-fr@enzolifesciences.com](mailto:info-fr@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)