

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 α 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 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

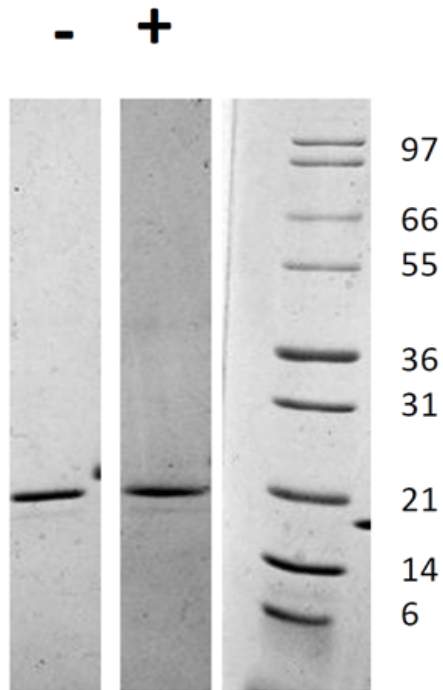


Figure 1: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human IL-6 has a predicted MW of 20.9 kDa.

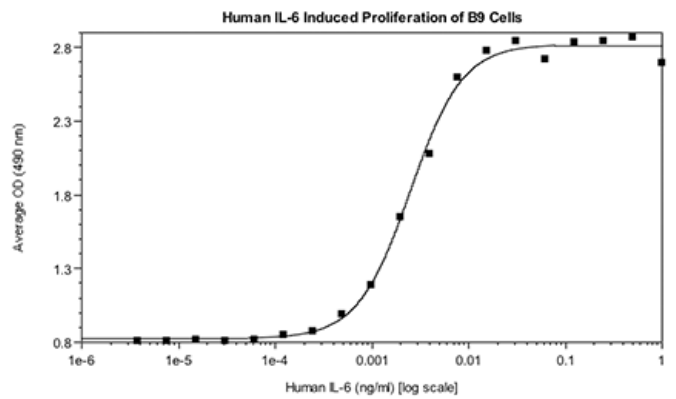


Figure 2: B9 cells were cultured with dilutions of Human IL-6. Cell proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

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- β 2, 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/ μ g 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	$\geq 97\%$ using Coomassie Staining
Reconstitution	Sterile water at 0.1 mg/mL
Source	Produced in <i>E. coli</i> . 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-
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)