

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 |
| ALX-201-818-0100 | 100 μ g |

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

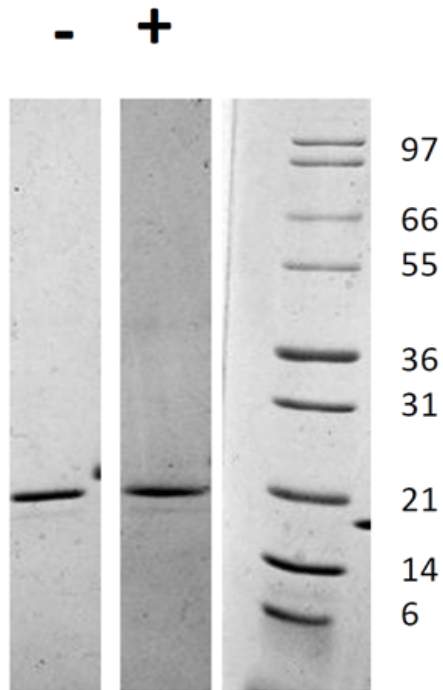


Figure 1: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human IL-6 has a predicted MW of 20.9 kDa.

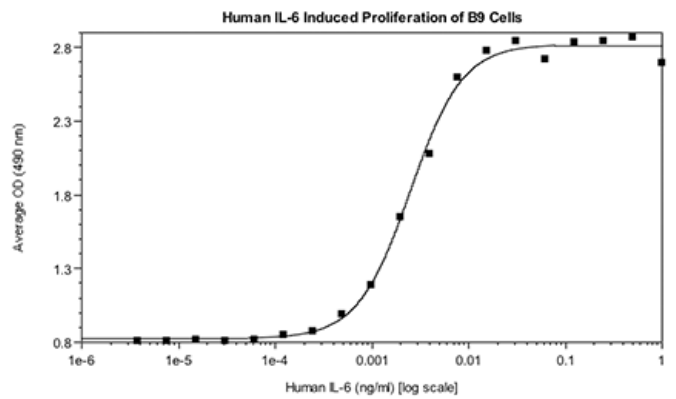


Figure 2: B9 cells were cultured with dilutions of Human IL-6. Cell proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

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-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 <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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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