# 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.

Carrier-free

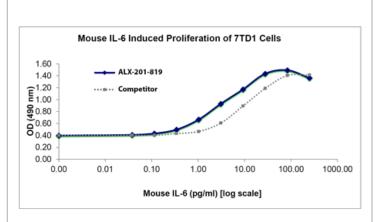
## **Ordering Information**

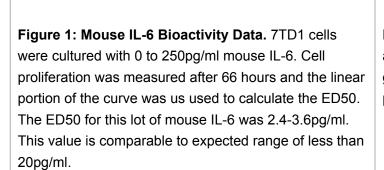
**Order Online** »

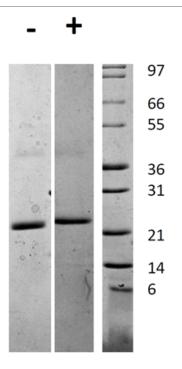
ALX-201-819-0010	10µg
ALX-201-819-0100	100µg

Manuals, SDS & CofA

**View Online** »







**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**

**Use/Stability** It is recommended that a carrier protein (0.1% HSA or BSA) is added for long term

storage.

**Handling** Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and

store at -20°C.

Long Term Storage -20°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 The activity is determined by dose-dependent induction of

7TD1 cell proliferation which is typically observed at

concentations < 0.02ng/ml.

Endotoxin Content ≤1 EU/µg protein measured by kinetic LAL analysis.

**Formulation** Lyophilized. Sterile filtered.

**MW** ~21.7kDa

**Purity** ≥95% (HPLC, Reducing and Non-reducing SDS-PAGE,

UV spectroscopy at 280 nm)

**Reconstitution** 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.

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

187 amino acids.

UniProt ID P08505

Last modified: May 29, 2024



info-

eu@enzolifesciences.com