

# Mouse IL-18 (recombinant)



**Alias:** IL-1F4 **Catalog #:** 6576

**Size:** 5 ug **Research Use Only**

**Molecular Weight:** 18.1 kDa

**Source:** Yeast. Recombinant Mouse IL-18 was produced in yeast and, therefore, does not have endotoxin. It is naturally folded and post-translationally modified.

**Formulation:** Lyophilized without carrier protein.

**Purity:** >95% as visualized by SDS-PAGE analysis.

**Purification:** Ion-exchange chromatography.

**Bioactivity:** In testing

**Entrez Gene ID:** 16173

**Number of Amino Acids:** 157

**Amino Acid Sequence:** NFGRLHCTTA VIRNINDQVL FVDKRQPVFE DMTDIDQSAS EPQTRLIIYM YKDSEVRGLA  
VTLSVKDSKM STLSCKNKII SFEEMDPPEN IDDIQSDLIF FQKRVPGHNK MEFESSLYEG  
HFLACQKEDD AFKLILKKKD ENGDKSVMFT LTNLHQS (157)

**Country of Origin:** USA

**Reconstitution:** Reconstitute with sterile phosphate-buffered saline containing at least 0.1% carrier protein.

**Stability and Storage:** Stable for up to twelve months from date of receipt at -20°C. Stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C. Avoid repeated freeze/thaw cycles.

**Applications:** The Mouse IL-18 protein can be used in cell culture, as an IL-18 ELISA Standard, and as a Western Blot Control.

**Background:** Interleukin-18 (IL-18) is a cytokine produced by macrophages and other cells that belongs to the IL-1 superfamily. IL-18 works together with IL-12 to induce cell-mediated immunity following infection with microbial products like lipopolysaccharide (LPS). After stimulation with IL-18, natural killer (NK) cells and certain T cells release another important cytokine called interferon-gamma (IFN-gamma) or type II interferon that plays an important role in activating the macrophages or other cells. Apart from its physiological role, IL-18 is also able to induce severe inflammatory reactions, which suggests its role in certain inflammatory disorders. The protein encoded by this gene is a proinflammatory cytokine. This cytokine can induce the IFN-gamma production of T cells. The combination of this cytokine and IL-12 has been shown to inhibit IL-4 dependent IgE and IgG1 production, and enhance IgG2a production of B cells.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse IL-18 updated on: 1/29/2014