

# Mouse IL-10 (recombinant)



**Alias:** none **Catalog #:** 6503

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

**Molecular Weight:** 18.8 kDa

**Source:** Yeast. Recombinant Mouse IL-10 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:** 16153

**Number of Amino Acids:** 160

**Amino Acid Sequence:**

SRGQYSREDNNCTHFPVGGQSHMLLELRTAFSQQVTKFFQTKDQLDNILLTDSLMLQDFKGYLGCCQALSEMI  
QFYLVEVMPQAEKHGPEIKEHLNSLGEKLTLMRLRRCHRFLPCENKSKAVEQVKSDFNKLQDQGV  
YKAMNEFDIFINCIEAYMMIKMKS (160)

**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-10 protein can be used in cell culture, as a IL-10 ELISA Standard, and as a Western Blot Control.

**Background:** The IL-10 family of cytokines consists of nine members: IL-10, IL-19, IL-20, IL-22, IL-24, IL-26, IL-28A, IL-28B, and IL-29. These cytokines elicit diverse host defense mechanisms. IL-10 family cytokines are essential for maintaining the integrity and homeostasis of tissue epithelial layers. By promoting innate immune response, members of this family can limit the damage caused by viral and bacterial infections. They can also facilitate the tissue-healing process in injuries caused by infection or inflammation. IL-10 itself is an anti-inflammatory cytokine. IL-10 family cytokines have indispensable functions in many infectious and inflammatory diseases.

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