

# Mouse CCL3 (recombinant)



**Alias:** MIP-1 alpha **Catalog #:** 6655

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

**Molecular Weight:** 7.9 kDa

**Source:** Yeast. Recombinant Mouse CCL3 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:** 20302

**Number of Amino Acids:** 69

**Amino Acid Sequence:** APYGADTPTA CCFSYSRKIP RQFIVDYFET SSLCSQPGVI FLTKRNRQIC ADSKETWVQE  
YITDLELNA (69)

**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 CCL3 protein can be used in cell culture, as a CCL3 ELISA Standard, and as a Western Blot Control.

**Background:** Chemokine ligand 3 (CCL3), also known as Macrophage inflammatory protein-1 alpha (MIP-1 alpha) is a small cytokine belonging to the CC chemokine family. There are at least 27 distinct members of the C-C subgroup reported for mammals. They are characterized by two adjacent cysteines. CC chemokines induce the migration of monocytes and other cell types such as NK cells and dendritic cells. CCL3/MIP-1 alpha is a chemoattractant for several different leukocytes, with varying degrees of potency (monocytes = T-cells > neutrophils > eosinophils).

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse CCL3 updated on: 6/23/2015