

Rat CCL4 (recombinant)



Alias: MIP-1 beta **Catalog #:** 6567

Size: 5 ug **Research Use Only**

Molecular Weight: 7.8 kDa

Source: Yeast. Recombinant Rat CCL4 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: 116637

Number of Amino Acids: 69

Amino Acid Sequence: APIGSDPPTS CCFSYTSRKI HRNFVMDYYE TSSLCSQPAV VFLTKKGRQI CADPSEPWVN
EYVNDLELN (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 Rat CCL4/MIP-1 beta protein can be used in cell culture, as an CCL4/MIP-1 beta ELISA Standard, and as a Western Blot Control.

Background: Chemokine ligand 4 (CCL4) is a small cytokine belonging to the CC chemokine family that is commonly known as MIP-1 beta. 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. CCL4 (MIP-1 beta) is a chemoattractant for natural killer cells, monocytes and a variety of other immune cells. CCL4 is involved in several inflammatory and autoimmune diseases including viral infection such as HIV-1/AIDS.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Rat CCL4 updated on: 1/29/2014