

Bovine CCL2 (recombinant)



Alias: MCP-1 **Catalog #:** 6383

Size: 5 ug **Research Use Only**

Molecular Weight: 8.8 kDa

Source: Yeast. Recombinant Bovine CCL2 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: 281043

Number of Amino Acids: 76

Amino Acid Sequence: QPDAINSQVA CCYTFNSKKI SMQRLMNYRR VTSSKCPKEA VIFKTILGKE LCADPKQKWV
QDSINYLNKK NQTPKP

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 Bovine MCP-1 protein can be used in cell culture, as an MCP-1 ELISA Standard, and as a Western Blot Control.

Background: Chemokine (C-C motif) ligand 2 (CCL2), also known as monocyte chemoattractant protein-1 (MCP-1), 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. CCL2 (MCP-1) recruits monocytes, memory T cells, and dendritic cells to sites of tissue injury and infection. CCL2 (MCP-1) is implicated in pathogenesis of several diseases characterized by monocyte infiltration, such as psoriasis, rheumatoid arthritis and atherosclerosis.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Bovine CCL2 updated on: 1/29/2014