

Bovine CXCL10 (recombinant)



Alias: IP-10 **Catalog #:** 6409

Size: 5 µg **Research Use Only**

Molecular Weight: 9.3 kDa

Source: Yeast. Recombinant Bovine CXCL10 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: 615107

Number of Amino Acids: 83

Amino Acid Sequence:

VPLSRNTRCSCIEISNGSVNPRSLEKLEVIPASQSCPRVEIIATMKKNGEKRLNPESKTIKNULLKAINK
QRTKRSPRTRKEA

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 IP-10/CXCL10 protein can be used in cell culture, as an IP-10/CXCL10 ELISA Standard, and as a Western Blot Control.

Background: C-X-C motif chemokine 10 (CXCL10) also known as Interferon gamma-induced protein 10 (IP-10) or small-inducible cytokine B10 is a member of the C-X-C chemokine family. CXCL10 (IP-10) is secreted by several cell types in response to IFN-gamma. These cell types include monocytes, endothelial cells and fibroblasts. CXCL10 (IP-10) has been attributed to several roles, such as chemoattraction for monocytes/macrophages, T cells, NK cells, and dendritic cells, promotion of T cell adhesion to endothelial cells, antitumor activity, and inhibition of bone marrow colony formation and angiogenesis. There have been 17 different C-X-C chemokines described in mammals, that are subdivided into two categories: those with a specific amino acid sequence (or motif) of glutamic acid-leucine-arginine (or ELR for short) immediately before the first cysteine of the C-X-C motif (ELR-positive), and those without an ELR motif (ELR-negative). ELR-positive C-X-C chemokines specifically induce the migration of neutrophils, and interact with chemokine receptors CXCR1 and CXCR2. C-X-C chemokines that lack the ELR motif are chemoattractant for lymphocytes. CXCL10 (IP-10) elicits its effects by binding to the cell surface chemokine receptor CXCR3.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Bovine CXCL10 updated on: 1/29/2014