

Mouse CXCL12 (recombinant)



Alias: SDF-1 **Catalog #:** 6508

Size: 5 ug **Research Use Only**

Molecular Weight: 11.6 kDa

Source: Yeast. Recombinant Mouse CXCL12 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: 20315

Number of Amino Acids: 99

Amino Acid Sequence: GKPVLSYRC PCRFFESHIA RANVKHLKIL NTPNCALQIV ARLKNNNRQV CIDPKLKWIQ
EYLEKALNKG RREEKVGKKE KIGKKKRQKK RKAQKRKN (99)

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 CXCL12 protein can be used in cell culture, as a CXCL12 ELISA Standard, and as a Western Blot Control.

Background: CXCL12 (formerly stromal cell-derived factor 1) was originally described as a vital chemoattractant for B cells and monocytes. CXCL12 is unlike other chemokines in that it binds to one single chemokine receptor, CXCR4, which itself is recognized by no other chemokines. CXCL12 binds to its receptor CXCR4 to mediate cell-type specific physiological processes including cellular migration, survival, and apoptosis. CXCL12 is produced physiologically in various tissues, and its receptor CXCR4 is also expressed on various haematopoietic and non-haematopoietic cells. CXCL12 and CXCR4 are essential for life as mice deficient in either gene are unable to survive much past birth. CXCL12 and CXCR4 play a role in many different diseases, including cancer, HIV, and rheumatoid arthritis.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse CXCL12 updated on: 5/23/2016