

Human CXCL9 (recombinant)



Alias: MIG **Catalog #:** 6565

Size: 5 µg **Research Use Only**

Molecular Weight: 11.7 kDa

Source: Yeast. Recombinant Human CXCL9 (MIG) 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: 4283

Number of Amino Acids: 103

Amino Acid Sequence: TPVVRKGRCS CISTNQGTH LQSLKDLKQF APSPSCEKIE IIATLKNGVQ TCLNPDSADV
KELIKKWEKQ VSQKKKQKNG KKHQKKKVLK VRKSQRSRQK KTT (103)

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 human CXCL9/MIG protein can be used in cell culture, as a CXCL9/MIG ELISA Standard, and as a Western Blot Control.

Background: Chemokine (C-X-C motif) ligand 9 (CXCL9), also commonly known as monokine induced by interferon-gamma (MIG), is a small cytokine belonging to the CXC chemokine family and is closely related to two other ELR-negative CXC chemokines: CXCL10 and CXCL11.

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.

CXCL9 (MIG) is a T-cell chemoattractant. Induced by IFN-gamma (IFN-γ), the ELR-negative chemokine CXCL9 (MIG) 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 Human CXCL9 (MIG) updated on: 2/5/2014