

Recombinant Human CXCL12 / SDF-1 α

CXCL12 is a small cytokine protein, which during embryo development, directs the migration of hematopoietic cells from fetal liver to bone marrow and also directs the formation of large blood vessels. In adults, CXCL12 plays an important role in angiogenesis by recruiting endothelial progenitor cells from the bone marrow through a CXCR4 dependent mechanism.

Materials Provided

Catalog Number:	92-1011
Quantity Supplied:	1 vial(s), 2 μ g per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	8 kg/mol
Source:	E. coli
Purity:	\geq 98% by SDS-PAGE and HPLC
Endotoxin Level:	< 0.1 ng/ μ g (<1EU/ μ g).
Formulation:	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
Storage:	\leq -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	20 μ M
Reconstitution Volume:	0.013 mL
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com