

Recombinant Murine CXCL16

Murine CXCL16 is a cytokine protein expressed by macrophages, monocytes, B cells, and Schwann cells, and expression is induced by IFNG and TNF- α . CXCL16 is also an integral protein that gets cleaved by ADAM10 to release the soluble active polypeptide into the extracellular space, and the soluble protein is chemotactic for activated CD8 T cells, inducing the process of calcium flux and regulating the migration of T cells. CXCL16 is an agonist for the CXCR6 receptor, but also serves as a receptor for oxidized LDL and phosphatidylserine.

Materials Provided

Catalog Number:	92-1021
Quantity Supplied:	1 vial(s), 5 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:	9900 g/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:	0.025 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