

Recombinant Murine CXCL13

Murine CXCL13 chemoattracts B cells and is one of the key players for B cell entry into lymphoid follicles. It takes part in the formation of ectopic lymphoid follicles. Murine CXCL13 chemoattracts B cells and is one of the key cytokines for B cell entry to lymphoid follicles. It takes part in the formation of ectopic lymphoid tissue in chronic inflammation, and elicits its functions primarily by interacting with its chemokine receptor, CXCR5.

Materials Provided

| | |
|-------------------------------|--------------------------|
| Catalog Number: | 92-1020 |
| Quantity Supplied: | 1 vial(s), 5 µg 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: | 9800 g/mol |
| Source: | E. coli |
| Purity: | ≥ 98% by SDS-PAGE and HPLC |
| Endotoxin Level: | < 0.1 ng/µg (<1EU/µg). |
| Formulation: | Sterile filtered through a 0.2 micron filter. Lyophilized with no additives |
| Storage: | ≤ -20°C Please avoid multiple freeze/thaw cycles. |

Reagent Preparation

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

| | |
|--------------------------------|-----------------------|
| Stock Concentration: | 25 µM |
| Reconstitution Volume: | 0.020 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