

Recombinant Murine CCL19

CCL19 chemoattracts naïve T cells and maturing dendritic cells into the T zones of secondary lymphoid organs. CCL19 controls the cytoarchitecture, rate of endocytosis, survival, apoptosis, migratory speed, and maturation of dendritic cells by binding to this chemokine receptor, CCR7.

Materials Provided

| | |
|-------------------------------|--------------------------|
| Catalog Number: | 92-1035 |
| 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: | 9200 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: | 10 µM |
| Reconstitution Volume: | 0.054 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