

Recombinant Human CCL20

CCL20 is a cytokine protein involved in the formation and function of mucosal lymphoid tissues via chemoattraction of lymphocytes and dendritic cells towards the epithelial cells surrounding these tissues. CCL20 also carries out its effects on its target cells by binding and activating the chemokine receptor, CCR6.

Materials Provided

Catalog Number:	92-1005
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:	8 kg/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:	20 µM
Reconstitution Volume:	0.021 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