

## Recombinant Murine MIP-3 $\alpha$ (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

<b>Catalog Number:</b>	<b>92-1204</b>
<b>Quantity Supplied:</b>	1 vial(s), 5 $\mu$ g per vial
<b>Reconstitution Buffer:</b>	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

<b>Molecular Weight:</b>	7900 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	$\geq$ 98% by HPLC
<b>Endotoxin Level:</b>	<0.01 EU per 1 $\mu$ g
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	10 $\mu$ M
<b>Reconstitution Volume:</b>	0.063 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)