

## Recombinant Murine CCL3

Murine CCL3 is a small cytokine protein that imparts a critical role in the recruitment of leukocytes to the site of inflammation. While both CCL3 and CCL4 attracts monocytes, macrophages, and dendritic cells to the site of infection, CCL3 attracts CD8+ T cells, while CD4+ T cells are more responsive to CCL4. In addition to these functions, CCL3 also induces inflammatory cytokine secretion, mast cell degranulation, and NK cell activation.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1015</b>
<b>Quantity Supplied:</b>	1 vial(s), 2 µg 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>	7800 g/mol
<b>Source:</b>	E. coli
<b>Purity:</b>	≥ 98% by SDS-PAGE and HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/µg (<1EU/µg).
<b>Formulation:</b>	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
<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 µM
<b>Reconstitution Volume:</b>	0.026 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)