

Recombinant Human CCL3

CCL3 is a cytokine protein involved in the acute inflammatory response for the recruitment and activation of polymorphonuclear leukocytes.

Materials Provided

Catalog Number:	92-1002
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:	7800 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:	20 µM
Reconstitution Volume:	0.032 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