

Recombinant Murine CCL1/I-309/TCA-3

CCL1 is a cytokine protein secreted by activated T cells to attract monocytes, NK cells, and immature B cells by binding with CCR8 cell surface receptors. It is considered to be one of the many chemokines having its genes located on chromosome 17.

Materials Provided

Catalog Number:	92-1199
Quantity Supplied:	1 vial(s), 25 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:	Recombinant
Purity:	> 95% by SDS-PAGE
Endotoxin Level:	<0.01 EU per 1µg
Formulation:	N/A
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:	0.160 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