

Recombinant Murine CCL25/TECK

CCL25 is a cytokine protein that is believed to play a role in the development of T cells. CCL25 is a chemotactic for thymocytes, macrophages, and dendritic cells, and elicits its response by binding to the chemokine receptor, CCR9.

Materials Provided

Catalog Number:	92-1207
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:	14 kg/mol
Source:	Recombinant
Purity:	≥ 97% by SDS-PAGE
Endotoxin Level:	<0.10EU 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:	0.089 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