

Recombinant Human CCL13

CCL13 is a cytokine protein involved in the induction of chemotaxis in monocytes, eosinophils, T lymphocytes, and basophils by binding to cell surface receptors like CCR2, CCR3, and CCR5. Activity of this protein has been linked to allergic reactions such as asthma, and can be induced by the inflammatory cytokines, including interleukin-1 and TNF- α .

Materials Provided

Catalog Number:	92-1003
Quantity Supplied:	1 vial(s), 5 μ g 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:	8600 g/mol
Source:	E. coli
Purity:	\geq 98% by SDS-PAGE and HPLC
Endotoxin Level:	< 0.1 ng/ μ g (<1EU/ μ g).
Formulation:	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
Storage:	\leq -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	25 μ M
Reconstitution Volume:	0.023 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