

Recombinant Human MCP-4 (CCL13)

Recombinant human protein expressed in E. coli cells

Catalog # M824-40N

Lot # G3211-91

Product Description

Recombinant Human MCP-4 (CCL13) was expressed in E. coli cells. The protein accession number is [Q99616](#)

Gene Aliases

Monocyte chemoattractant protein 4, CCL13, NCC-1, C-C motif chemokine 13, Small-inducible cytokine A13, MCP4, NCC1, SCYA13, Recombinant Human Monocyte Chemoattractant Protein-4 (CCL13)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

Formulation

Lyophilized from a 0.2 µm filtered solution in PB, 100mM NaCl, pH 7.5

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

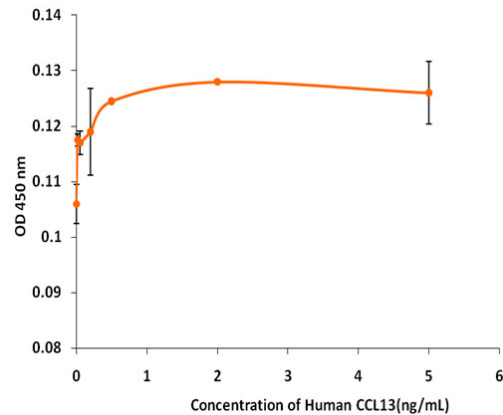
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

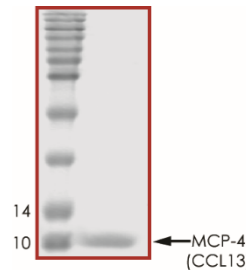
Monocyte Chemoattractant Proteins 4 (MCP-4/CCL13) is member of a distinct, structurally-related subclass of CC chemokines. MCP-4/CCL13 is a major chemoattractants for eosinophils, basophils, monocytes, and T lymphocytes. The MCP protein family bind to specific G-protein-coupled receptors, initiating a signal cascade within the cell. Expression of MCP-4/CCL13 mRNA and protein is greater in the sputum, epithelium, submucosal inflammatory cells, and bronchoalveolar lavage fluid of asthmatics than in healthy individuals. Involvement of MCPs has also been demonstrated in renal inflammation and atopic dermatitis. The expression of this protein is also correlated with enhanced inflammatory immune responses during immunotherapy. The recombinant construct corresponds to the "short-chain" of CCL13 spanning amino acids 24 – 98.

Specific Activity



Determined by its ability to chemoattract human monocytes using a concentration range of 20-40 ng/mL.

Purity



Recombinant Human MCP-4 (CCL13) resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **9.0 kDa**

Recombinant Human MCP-4 (CCL13)

Recombinant human protein expressed in E. coli cells

Catalog Number	M824-40N
Specific Lot Number	G3211-91
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.