

## Recombinant Human MIP-4 (CCL18)

Recombinant human protein expressed in E. coli cells

**Catalog # M844-40N**

Lot # G3211-97

### Product Description

Recombinant Human MIP-4 (CCL18) was expressed in E. coli cells. The protein accession number is [P55774](#)

### Gene Aliases

MIP4, CCL18, Pulmonary and activation-regulated chemokine, PARC, DC-CK1, AMAC-1, C-C motif chemokine 18, Small-inducible cytokine A18, Recombinant Human Macrophage Inflammatory Protein-4 (CCL18)

### Endotoxin Level

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

### Formulation

Recombinant CC18 was lyophilized from a 0.2 µm filtered 20 mM PB, 100 mM NaCl solution 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.

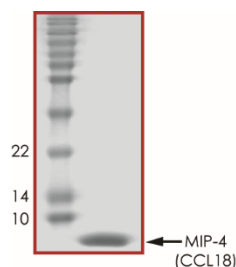
### Scientific Background

CCL18, also designated dendritic cell (DC)3-derived CC chemokine 1, pulmonary and activation-regulated chemokine, alternative macrophage activation-associated CC chemokine 1, and MIP-4, is a human chemokine structurally related to CCL3. CCL18 was shown to be expressed in germinal centers of tonsils by dendritic cells and to attract mainly naive T cells, CD38 mantle zone B lymphocytes and DC. The production of CCL18 by APC is enhanced by Th2 cytokines like IL-4 and IL-13 and suppressed by IFN-gamma.

### Specific Activity

Determined by its ability to chemoattract human T cells using a concentration range of 2-40 ng/mL.

### Purity



Recombinant Human MIP-4 (CCL18) resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **8.0 kDa**

## Recombinant Human MIP-4 (CCL18)

Recombinant human protein expressed in E. coli cells

Catalog Number	M844-40N
Specific Lot Number	G3211-97
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.
	Usage

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

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