



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

MIP-3 alpha (CCL20)

Catalog #: 90220-B (25 µg)
Lot #: 150723

Description: Macrophage inflammatory protein (MIP)-3/CCL20, also known as liver and activation-regulated chemokine (LARC) or Exodus, is a member of the CC chemokine subfamily initially noted to be expressed in human liver, lung, appendix, and tonsillar crypts. MIP-3 is selectively chemotactic for CD34+ bone marrow cell-derived immature DCs and CD45RO+ memory T cells that express the cognate receptor CCR6. MIP-3 produced at sites of inflammation may chemoattract CCR6-expressing immature DCs to the subepithelial region of mucosal surfaces.

Source: Optimized DNA sequence encoding human MIP-3 alpha mature chain was expressed in *E. coli*.

Formulation: Lyophilized from a 0.2 µm filtered solution containing 20 mM phosphate buffer, pH 7.5, 100 mM NaCl.

Reconstitution: Reconstitute at 0.1-1.0 mg/ml in distilled water. This solution can then be diluted into other buffers. To maximize product collection from vial surface, vortex briefly and then spin down to recollect the liquid.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >98%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/µg (1EU/µg), using the LAL gel clot method.

Biological Activity: Determined by its ability to chemo-attract human monocytes using a concentration range of 2.0-40.0 ng/ml.

Protein Sequence:
ASNFDCCLGYTDRILHPKFIVGFTRQL
ANEGCDINAIIFHTKKKLSVCANPKQT
WVKYIVRLLSKKVKNM

References:

1. J Rheumatol, Nov 2009, 36: 2397 - 2402.2.
2. Arterioscler Thromb Vasc Biol, Oct 2009, 29: 1608 - 1614.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com