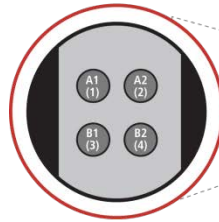


MSD[®] Mouse MIP-3 α Kit

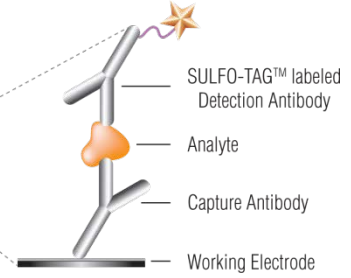
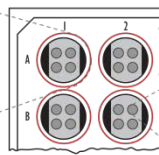
For quantitative determination in mouse serum and plasma



1. MIP-3 α
2. BSA blocked
3. BSA blocked
4. BSA blocked



MSD MULTI-SPOT[®]
96-Well 4-Spot Plate



Macrophage inflammatory protein 3 alpha (MIP-3 α) (LARC/CCL20) is a C-C chemokine with inflammatory and homeostatic functions.^{1,2} Initially identified in the liver, it is expressed in lymphatic tissue, lung tissue, macrophages, dendritic cells, B- and T- lymphocytes, and eosinophilic granulocytes as well as in normal colon, pancreas, prostate, uterine cervix, and skin.² To date, MIP-3 α is the only known ligand for the CCR6 receptor; it is also chemotactic for CCR6⁺ cells such as Th17.¹⁻³

MIP-3 α is implicated in a broad spectrum of disorders, including colorectal cancer and tumor metastasis,¹ rheumatoid arthritis,³ psoriasis,⁴ obesity,⁵ and wound healing.⁶ Disrupting the MIP-3 α and CCR6 interaction may ultimately prove to be a viable therapeutic strategy,^{4,7} as CCR6 deletion resulted in reduced atherosclerotic lesion area and reduced macrophage presence at atherosclerotic plaque sites in models of atherogenesis.⁷

The assay is available on 96-well, 4-spot plates. Representative data from the assay is presented below.

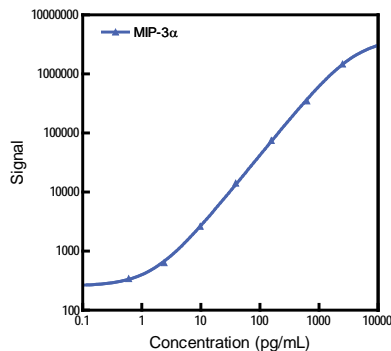
Alzheimer's Disease
BioProcess
Cardiac
Cell Signaling
Clinical Immunology
Cytokines
Growth Factors
Hypoxia
Immunogenicity
Inflammation
Metabolic
Oncology
Toxicology
Vascular

Catalog Numbers

Mouse MIP-3 α Kit	
Kit Size	Catalog #
1 plate	K152MSD-1
5 plates	K152MSD-2
25 plates	K152MSD-4

Assay Sensitivity

The following standard curve illustrates the dynamic range of the Mouse MIP-3 α assay.



MIP-3 α	
Average LLOD (pg/mL)	0.33

The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator blank).

Specificity

To assess specificity of the MIP-3 α assay, the kit was tested with the following recombinant mouse proteins: IFN γ , IL-1 β , IL-2, IL-4, IL-5, IL-6, KC, IL-10, IL-12, TNF α , IL-23, GM-CSF, IL-13, VEGF, RANTES, TNF-R1, TNF-R2, IL-12/IL-23 p40, and MCP-1 at 1250 pg/mL. Less than 0.1% non-specific binding was observed with each protein.

Ordering Information

MSD Customer Service
Phone: 1-301-947-2085
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850 USA
www.mesoscale.com[®]

For Research Use Only.
Not for use in diagnostic procedures.

MSD Cytokine Assays

MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25 μ L or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference
- **Simple protocols:** Only labels bound near the electrode surface are excited, enabling assays with fewer washes
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple rounds of label excitation and emission enhance light levels and improve sensitivity

For a complete list of products, please visit our website at www.mesoscale.com.

References

1. Schutyser E, et al. Regulated production and molecular diversity of human liver and activation-regulated chemokine/macrophage inflammatory protein-3 alpha from normal and transformed cells. *J Immunol.* 2000 Oct 15;165(8):4470-7.
2. Ghadjar P, et al. The chemokine CCL20 and its receptor CCR6 in human malignancy with focus on colorectal cancer. *Int J Cancer.* 2009 Aug 15;125(4):741-5.
3. Hirota K, et al. Preferential recruitment of CCR-6 expressing Th17 cells to inflamed joints via CCL20 in rheumatoid arthritis and its animal model. *J Exp Med.* 2007 Nov 26;204(12):2803-12.
4. Harper EG, et al. Th17 cytokines stimulate CCL20 expression in keratinocytes *in vitro* and *in vivo*: implications for psoriasis pathogenesis. *J Invest Dermatol.* 2009 Sep;129(9):2175-83.
5. Duffant C, et al. Interplay between human adipocytes and T lymphocytes in obesity: CCL20 as an adipochemokine and T lymphocytes as lipogenic modulators. *Arterioscler, Thromb, Vasc Biol.* 2009 Oct;29(10):1608-14.
6. Kennedy-Crispin M, et al. Human keratinocytes' response to injury upregulates CCL20 and other genes linking innate and adaptive immunity. *J Invest Dermatol.* 2012 Jan;132(1):105-13.
7. Wan W, et al. Genetic deletion of chemokine receptor Ccr6 decreases atherogenesis in ApoE-deficient mice. *Circ Res.* 2011 Aug 5;109(4):374-81.
8. Hem A, et al. Saphenous vein puncture for blood sampling of the mouse, rat, hamster, gerbil, ferret and mink. *Lab Anim.* 1998 Oct;32(4):364-8.
9. Removal of blood from laboratory animals and birds: First report of the BVA/FRAME/RSPCA/UFAW joint working group on refinement. *Lab Animal* 1993 Jan;27(1):1-22.

