

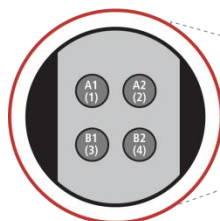
MSD® Human MIP-4 Kit

For quantitative determination in human serum and plasma

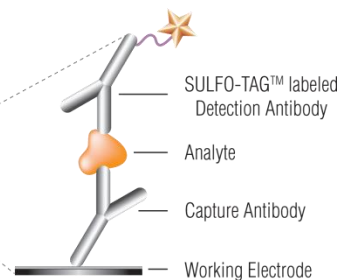
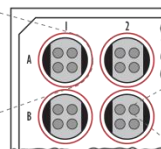


Alzheimer's Disease
BioProcess
Cardiac
Cell Signaling
Clinical Immunology
Cytokines
Growth Factors
Hypoxia
Immunogenicity
Inflammation
Metabolic
Oncology
Toxicology
Vascular

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



MSD MULTI-SPOT®
96-Well 4-Spot Plate



Macrophage inflammatory protein 4 (MIP-4) (CCL18/PARC/AMAC-1/DC-CK-1/SCYA18) is a C-C chemokine strongly expressed in the lung and to a lesser extent in the thymus and lymph nodes.^{1,2} To date, the functional receptor and role of MIP-4 has been difficult to ascertain as it appears to exist only in primates, rendering rodent models unusable. MIP-1 α is the most closely related chemokine to MIP-4, but CCR1 and CCR5 (receptors for MIP-1 α) do not bind MIP-4.¹ *In vitro*, MIP-4 is particularly chemotactic for naive T-lymphocytes, Th2 cells, B cells, and immature dendritic cells.^{1,2} Despite MIP-4's absence in rodents, it is functionally active in humans as a T-lymphocyte chemoattractant.³

Research implicates MIP-4 involvement in tumor malignancy and various inflammatory and allergic pulmonary, skin, and joint diseases.⁴ In addition to its chemotactic function, MIP-4 is able to induce production of adaptive regulatory T cells from CD4⁺CD25⁺ T cells.² However, this phenomenon is only present in individuals with non-allergic asthma. Taken in conjunction with elevated MIP-4 levels in allergic diseases (asthma and atopic dermatitis), this suggests MIP-4 desensitization or decreased feedback regulation.² In chronic obstructive pulmonary disease, an elevated level of MIP-4 coincides with clinical outcome and can serve as a potential biomarker.⁵ MIP-4 is also expressed in abundance in tumor-associated macrophages found in breast tissue and is associated with increased breast cancer metastasis.⁶ MIP-4 is thought to cause increased ECM adherence of cancerous cells to fibronectin in the stroma allowing subsequent invasion.⁵

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

Catalog Numbers

Human MIP-4 Kit	
Kit Size	Catalog #
1 plate	K151RLD-1
5 plates	K151RLD-2
25 plates	K151RLD-4

Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

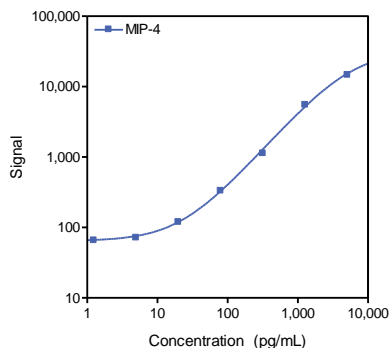
Company Address

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.

Assay Sensitivity

The following standard curve illustrates the dynamic range of the Human MIP-4 assay.



MIP-4	
Average LLOD (pg/mL)	9.06

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-4 assay, the kit was tested with the following recombinant human proteins: fractalkine, 35,000 pg/mL; I-TAC, 1,500 pg/mL; MCP-2, 250 pg/mL; MIP-3 β , 275 pg/mL; and MIP-5, 1,200 pg/mL. Less than 0.1% non-specific binding was observed with each protein.

MSD Cytokine Assays

MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25 µL 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. Schraufstatter IU, et al. The chemokine CCL18 causes maturation of cultured monocytes to macrophages in the M2 spectrum. *Immunology*. 2011 April;135(4):287-98.
2. Chang Y, et al. The chemokine CCL18 generates adaptive regulatory T cells from memory CD4⁺ T cells of healthy but not allergic subjects. *FASEB J*. 2010 Dec;24(12):5063-72.
3. Luzina IG; Atamas, SP. CCR6 is not necessary for functional effects of human CCL18 in a mouse model. *Fibrogenesis Tissue Repair*. 2012 Jan 18;5(1):2.
4. Schutyser E, et al. Involvement of CC chemokine ligand 18 (CCL18) in normal and pathological processes. *J Leukoc Biol*. 2005 Jul;78(1):14-26.
5. Sin DD, et al. Serum PARC/CCL-18 concentrations and health outcomes in chronic obstructive pulmonary disease. *Am J Respir Crit Care Med*. 2011 May 1;183(9):1187-92.
6. Chen J, et al. CCL18 from tumor-associated macrophages promotes breast cancer metastasis via PITPNM3. *Cancer Cell*. 2011 Apr 12;19(4):541-55.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, S-PLEX, U-PLEX, V-PLEX, MSD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), S-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2012, 2017 Meso Scale Diagnostics, LLC. All rights reserved.

