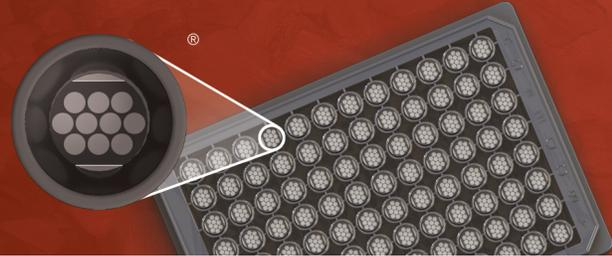


U-PLEX[®] Human MCP-2 Assay



www.mesoscale.com[®]

Ordering Information

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

Scientific Support

Phone: 1-301-947-2025
Email: ScientificSupport@mesoscale.com

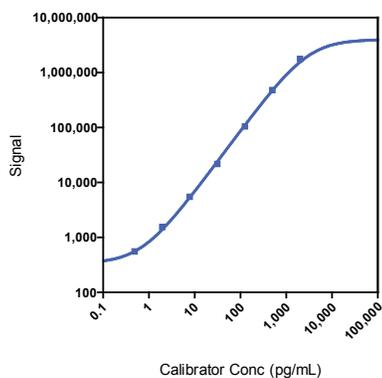
Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

| | |
|---------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Product Options | Available as part of U-PLEX Biomarker Group 1 (hu) multiplex combination: K15067L-1/-2/-4 |
| | Individual assay: K151XHK-1/-2/-4; Antibody Set: B21XH-2/B21XH-3 |
| | For more ordering options, please visit www.mesoscale.com |
| Instrument Compatibility | SECTOR [®] Imager 2400, SECTOR Imager 6000, MESO [®] SECTOR S 600, MESO QuickPlex [®] SQ 120 |
| Sample Type | Human serum, EDTA plasma, and cell culture supernatants |
| Assay Protocol | Refer to the U-PLEX Biomarker Group 1 (Human) product insert available at www.mesoscale.com/U-PLEX-documents |

The U-PLEX platform was designed to provide ultimate flexibility for detection of biomarkers in a wide variety of sample types. This datasheet provides the representative performance of the U-PLEX Human MCP-2 Assay tested on U-PLEX plates run as a multiplex. The data were generated during the development of the assay and do not represent the product specifications. Under your experimental conditions and with your specific multiplex, the assay may perform differently than the representative data shown. U-PLEX assays are available in multiplex format with other compatible assays. The same assay can also be used to detect a single analyte on MSD GOLD[™] Small Spot Streptavidin plates.

Representative Calibration Curve and Sensitivity



| Assay | Median LLOD (pg/mL) | LLOD Range (pg/mL) |
|-------|---------------------|--------------------|
| MCP-2 | 0.11 | 0.11-0.17 |

The calibration curves used to calculate analyte concentrations were established by fitting the signals from the Calibrators using a 4-parameter logistic (or sigmoidal dose-response) model with a $1/Y^2$ weighting. Analyte concentrations were determined from the electrochemiluminescence signals by back-fitting to the calibration curve. The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero Calibrator).

Precision

| | Control | Average Conc. (pg/mL) | Average Intra-run Conc. %CV | Inter-run Conc. %CV |
|-------|---------|-----------------------|-----------------------------|---------------------|
| MCP-2 | High | 251 | 4.0 | 10.4 |
| | Mid | 50.4 | 4.7 | 10.1 |
| | Low | 11.3 | 7.0 | 12.8 |

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

Controls were made by spiking Calibrator into assay diluent at 3 levels within the quantitative range of the assay. Average intra-run concentration %CV is the average %CV of the control replicates within an individual run. Inter-run concentration %CV is the variability of controls across multiple runs.

Note: This datasheet contains representative assay performance data. In custom multiplex formats, the assay may perform differently than the representative data shown.

MSD® U-PLEX Assays

Spike Recovery

| | Spike Level | Serum | | Plasma | | Cell Culture Media | |
|-------|-------------|--------------------|------------------|--------------------|------------------|--------------------|------------------|
| | | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range |
| MCP-2 | High | 85 | 75-97 | 82 | 56-94 | 104 | 98-108 |
| | Mid | 86 | 78-92 | 92 | 87-96 | 96 | 93-101 |
| | Low | 92 | 88-101 | 97 | 96-98 | 93 | 90-95 |

Normal human serum, EDTA plasma, and cell culture media were spiked with Calibrator at 3 levels. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may require additional dilution with assay diluent to reduce matrix effects.

% Recovery = (measured concentration / expected concentration) x 100

Tested Samples

| Sample Type | Serum | Plasma | Stimulated Cell Models |
|----------------|-------|--------|------------------------|
| Median (pg/mL) | 28 | 21 | 1.6 |
| Range (pg/mL) | 20-37 | 18-74 | ND-AS |
| % Detected | 100 | 100 | 80 |

ND = non-detectable (< LLOD), AS = above standard 1

Normal serum and EDTA plasma samples were tested without dilution prior to the assay.

Dilution Linearity

| | Serum | | | Plasma | | | Cell Culture Media | | |
|-------|---------------|--------------------|------------------|---------------|--------------------|------------------|--------------------|--------------------|------------------|
| | Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range |
| MCP-2 | 2 | 96 | 92-104 | 2 | 98 | 97-99 | 2 | 94 | 87-98 |
| | 4 | 91 | 87-97 | 4 | 93 | 89-95 | 4 | 90 | 79-104 |
| | 8 | 81 | 77-87 | 8 | 88 | 83-90 | 8 | 79 | 69-92 |

Normal human serum, EDTA plasma, and cell culture media were spiked with recombinant Calibrator and tested at different dilutions. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may require additional dilution with assay diluent to reduce matrix effects.

% Recovery = (measured concentration / expected concentration) x 100

Specificity

To assess specificity, the MCP-2 Antibody Set was tested individually against a larger panel of recombinant human analytes for nonspecific binding (CTACK, ENA-78, Eotaxin, Eotaxin-2, Eotaxin-3, EPO, FLT3L, Fractalkine, G-CSF, GM-CSF, GRO- α , I-309, IFN- α 2a, IFN- β , IFN- γ , IL-1 α , IL-1 β , IL-1RA, IL-2, IL-2R α , IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12/IL-23p40, IL-12p70, IL-13, IL-15, IL-16, IL-17A, IL-17A/F, IL-17B, IL-17C, IL-17D, IL-17E/IL-25, IL-17F, IL-18, IL-21, IL-22, IL-23, IL-27, IL-29/IFN- λ 1, IL-31, IL-33, IP-10, I-TAC, MCP-1, MCP-2, MCP-3, MCP-4, M-CSF, MDC, MIF, MIP-1 α , MIP-1 β , MIP-3 α , MIP-3 β , MIP-5, SDF-1 α , TARC, TNF- α , TNF- β , TPO, TRAIL, TSLP, VEGF-A, and YKL-40). Nonspecific binding was less than 0.5%. MCP-2 detection antibody nonspecifically binds (2.8%) to Eotaxin.

% Nonspecificity = (nonspecific signal / specific signal) x 100

Diluent Compatibility

The data included in this document have been collected using Diluents 3 and 43. MSD offers a range of assay and antibody diluents for separate purchase. Depending on your assay needs, other diluents may be tested.

Assay Components

Calibrator: Human MCP-2 is included in Calibrator 10 blend. The full-length recombinant human protein expressed in E.coli is used.

Antibodies: The U-PLEX Human MCP-2 Assay uses mouse monoclonal antibody for capture and rabbit polyclonal antibody for detection.

Assay generation: A

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), R-PLEX (design), S-PLEX (design), T-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.

©2016, 2018 Meso Scale Diagnostics, LLC. All rights reserved.

