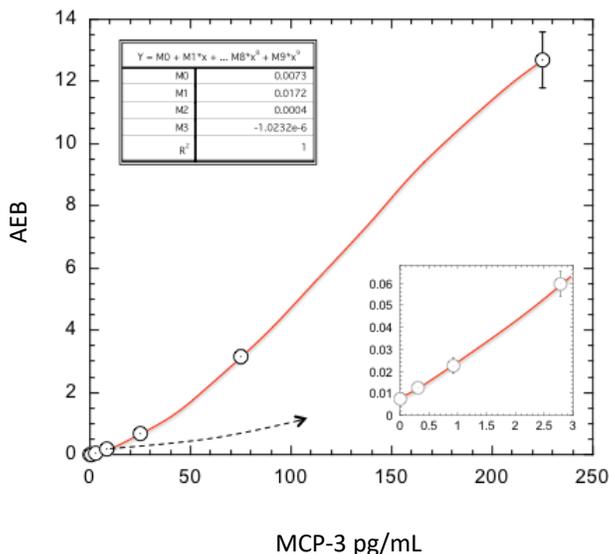


**Description**

Human monocyte chemotactic protein-3 is a 99 residue, 11 kDa secreted protein which shares 73% sequence similarity to its most structurally and functionally related homolog, MCP-1. MCP-3 displays strong interaction with CC chemokine receptor 2 (CCR2). Chemotactic gradients of MCP-3 are generated in response to localized cell infection which sequesters phagocytes to the site of inflammation and MCP-3 has been well implicated in the process of inflammation in response to presence of allergen. Elevated MCP-3 levels are associated with cytokine stimulated human osteosarcoma (cell line MG-63) and have been documented in patients with inflammatory bowel disease (IBD), ulcerative colitis, and Crohn’s disease. Faulty MCP-3 immune stimulation of non-inflamed lesioned areas of the CNS has been reported in patients with MS.

**Calibration Curve:** Cubic fit parameters are depicted.



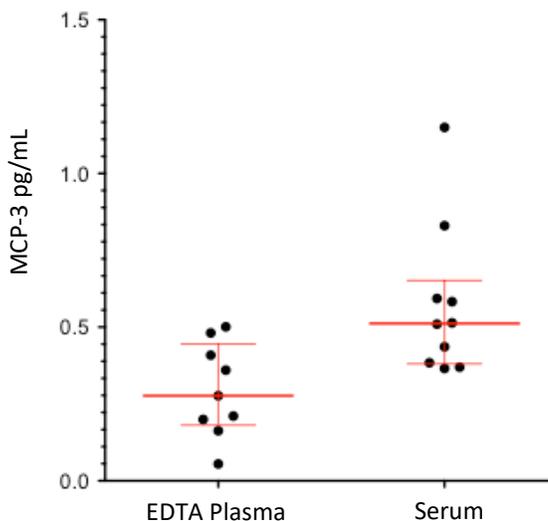
**Lower Limit of Quantification (LLOQ):** Triplicate measurements of serially diluted calibrator were read back on the calibration curve over 1 reagent lot on 1 instrument (5 runs total).

**Limit of Detection (LOD):** Calculated as 2.5 standard deviations from the mean of background signal read back on each calibration curve over 1 reagent lot on 1 instrument (5 runs total).

<b>LLOQ</b>	<b>0.309 pg/mL</b> pooled CV 15.9% mean recovery 104%
<b>LOD</b>	<b>0.124 pg/mL</b> range 0.0597–0.222 pg/mL
<b>Dynamic range (serum and plasma)</b>	0–450 pg/mL
<b>Diluted Sample volume*</b>	120 µL per measurement
<b>Tests per kit</b>	192

\*See Kit Instruction for details

**Endogenous Sample Reading:** Healthy donor matched EDTA plasma (n=9) and serum (n=10) were measured. 1 plasma sample was not detectable. Error bars depict median and interquartile ranges.



Sample Type	Median MCP-3 pg/mL	% Above LOD
EDTA Plasma	0.445	90%
Serum	0.512	100%

**Precision:** Representative precision was estimated with repeated assay of serum and plasma panels using two instruments and one reagent lot. Within-run and between-run CVs are depicted in the following table. Within-run CVs reflect average CVs across 5 experiments of 3 replicates each.

Sample	Mean (pg/mL)	Within run CV	Between run CV
Serum Panel 1	87.5	7.3%	5.1%
Serum Panel 2	326	3.9%	9.9%
Plasma Panel 3	6.91	10.2%	11.8%
Serum Panel 4	5.94	9.7%	9.0%
Plasma Panel 5	10.9	7.3%	8.7%

**Spike and Recovery:** MCP-3 spiked into 6 serum samples at 2 levels.

**Dilution Linearity:** Spiked, diluted 2x serially from MRD (2x) to 128x with Sample Diluent.

<b>Spike and Recovery (Serum)</b>	<b>Mean = 89.9%</b> Range: 46.7–146%
<b>Dilution Linearity (128x)</b>	<b>Mean = 118%</b> Range: 92.7–144%