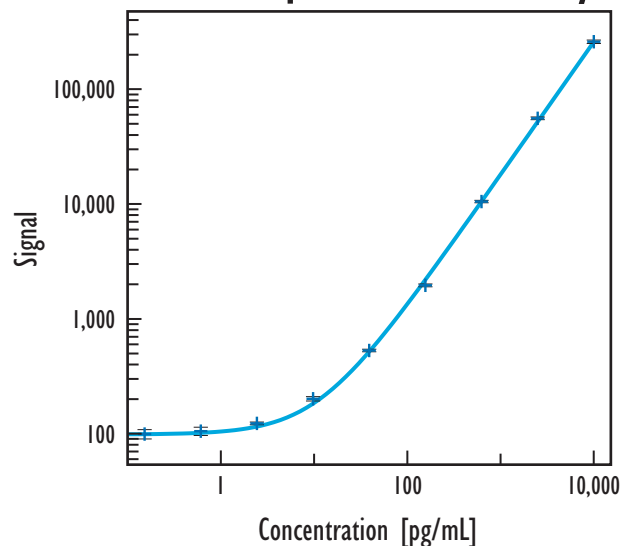


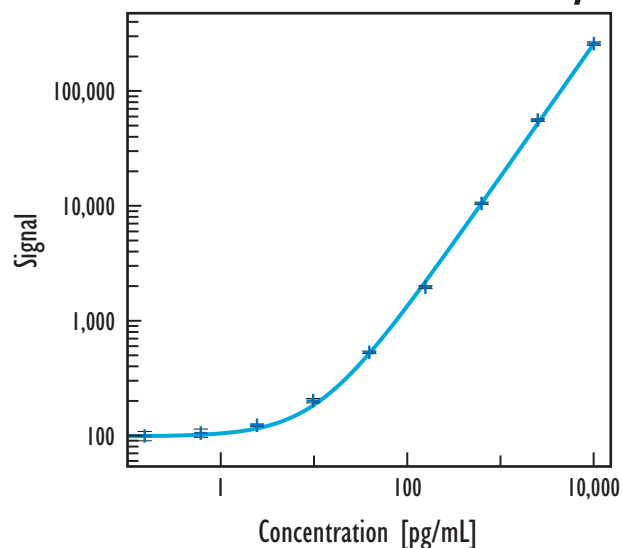
Mouse MCP-1 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse MCP-1		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	89	8.2
0.010	96	7.3
0.038	101	2.4
0.15	99	9.3
0.61	105	8.2
2.4	123	2.7
9.8	202	4.3
39	531	2.6
156	1,960	2.8
625	10,518	2.5
2,500	55,835	2.7
10,000	259,246	3.4
40,000	296,470	6.0

Serum / Plasma Assay



Mouse MCP-1		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	52	10.4
0.15	42	10.4
0.61	38	29.0
2.4	61	3.5
9.8	87	3.4
39	196	3.7
156	694	9.6
625	3,231	10.8
2,500	19,453	9.2
10,000	94,272	9.1
40,000	310,549	1.8

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	4.1
Serum / Plasma	9.0

Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	97
Serum	94
EDTA Plasma	97
Heparin Plasma	87

Spike recoveries were determined in each matrix over a range of spike levels from 39 to 625 pg/mL. Each spike was tested in triplicate.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	0	63 - 222
EDTA Plasma (n=2)	0	45 - 334
Heparin Plasma (n=10)	0	43 - 274

Endogenous cytokine levels (pg/mL) were determined for different sample types across multiple samples. (n/d indicates cytokine level below the detection limit)

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested.



Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.
www.mesoscale.com