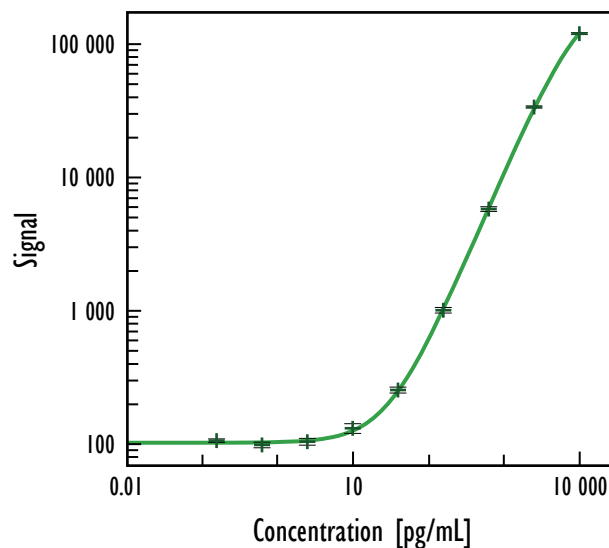


MULTI-ARRAY[®] Human MCP-1 Tissue Culture Assay

Detection of MCP-1 in Tissue Culture Samples

Typical Standard Curve Data



Human MCP-1	
Concentration (pg/mL)	Mean Signal
0	104
0.2	109
0.6	101
2	107
10	134
39	259
156	1 021
625	5 797
2 500	33 406
10 000	117 771

Average %CV	4.5
-------------	-----

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Tissue Culture	10

Definition of Detection Parameters

LLOD = Lower Limit of Detection is 2.5 stdev above the background

Standard curve data is from a representative experiment

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K151AYB-1
Tissue Culture	5 plates	K151AYB-2
Tissue Culture	25 plates	K151AYB-4
Ultra-Sensitive	1 plate	K151AYC-1
Ultra-Sensitive	5 plates	K151AYC-2
Ultra-Sensitive	25 plates	K151AYC-4

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

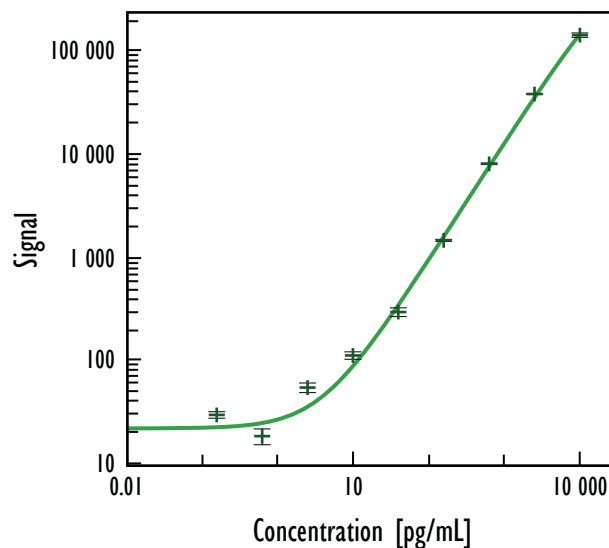


Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.
www.mesoscale.com

MULTI-ARRAY[®] Human MCP-1 Ultra-Sensitive Assay

Detection of MCP-1 in Human Serum and Plasma Samples

Typical Standard Curve Data



Human MCP-1	
Concentration (pg/mL)	Mean Signal
0	23
0.2	30
0.6	18
2	54
10	112
39	294
156	1 463
625	8 103
2 500	38 596
10 000	144 230
Average %CV	5.4

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Serum/Plasma	4.2

Definition of Detection Parameters

LLOD = Lower Limit of Detection is 2.5 stdev above the background

Standard curve data is from a representative experiment

Spike Recoveries

Sample Type	Average % Recovery
Serum	107
EDTA Plasma	116
Heparin Plasma	116

Spike recoveries were determined in each matrix over a range of spike levels.