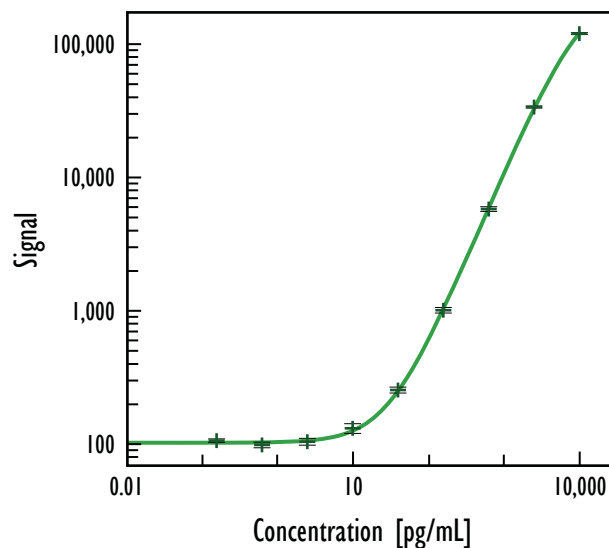


MULTI-ARRAY[®] Human MCP-1 Tissue Culture Assay

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

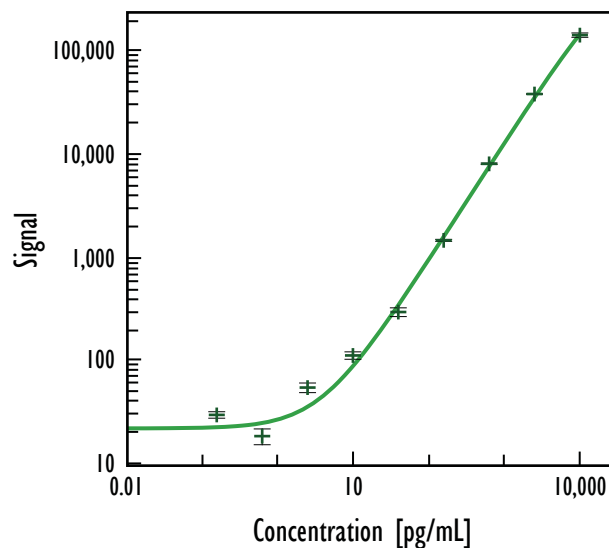
Standard curve data is from a representative experiment

Kit Model	Kit Size	SI2400	SI6000
Tissue Culture	1 plate	K151AYB-1	K111AYB-1
Tissue Culture	5 plates	K151AYB-2	K111AYB-2
Tissue Culture	25 plates	K151AYB-4	K111AYB-4
Ultra-Sensitive	1 plate	K151AYC-1	K111AYC-1
Ultra-Sensitive	5 plates	K151AYC-2	K111AYC-2
Ultra-Sensitive	25 plates	K151AYC-4	K111AYC-4
Base	25 plates	K151AYA-4	K111AYA-4
Base	250 plates	K151AYA-5	K111AYA-5

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.