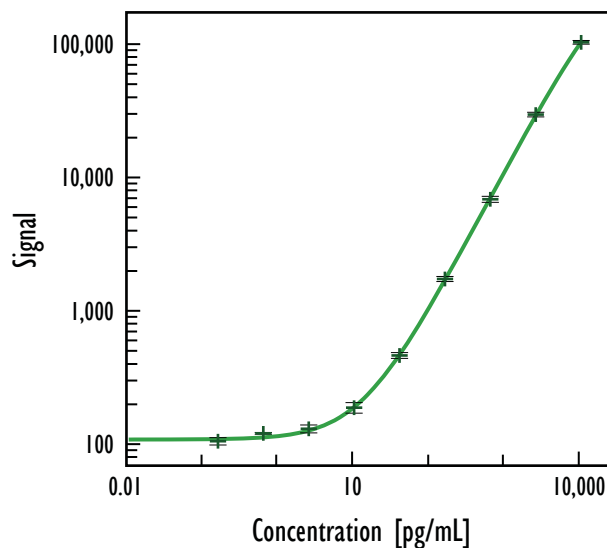


MULTI-ARRAY[®] Human MIP-1 β Tissue Culture Assay

Detection of MIP-1 β in Tissue Culture Samples

Typical Standard Curve Data



Human MIP-1 β	
Concentration (pg/mL)	Mean Signal
0	96
0.2	97
0.6	111
2	121
10	174
39	434
156	1,649
625	6,629
2,500	29,084
10,000	102,660

Average %CV	5.4
-------------	-----

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Tissue Culture	3.5

Definition of Detection Parameters

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

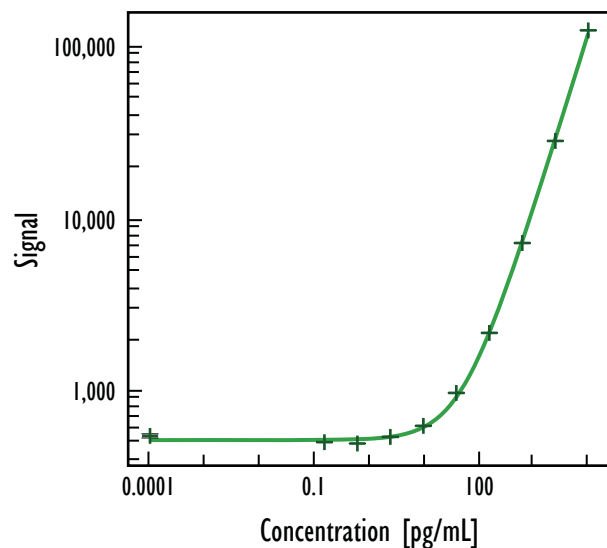
Standard curve data is from a representative experiment

Kit Model	Kit Size	SI2400	SI6000
Tissue Culture	1 plate	K151BDB-1	K111BDB-1
Tissue Culture	5 plates	K151BDB-2	K111BDB-2
Tissue Culture	25 plates	K151BDB-4	K111BDB-4
Ultra-Sensitive	1 plate	K151BDC-1	K111BDC-1
Ultra-Sensitive	5 plates	K151BDC-2	K111BDC-2
Ultra-Sensitive	25 plates	K151BDC-4	K111BDC-4
Base	25 plates	K151BDA-4	K111BDA-4
Base	250 plates	K151BDA-5	K111BDA-5

MULTI-ARRAY[®] Human MIP-1 β Ultra-Sensitive Assay

Detection of MIP-1 β in Human Serum and Plasma Samples

Typical Standard Curve Data



Human MIP-1 β	
Concentration (pg/mL)	Mean Signal
0	545
0.15	501
0.61	492
2.4	538
9.8	623
39	969
156	2,164
625	7,201
2,500	28,257
10,000	124,201

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Serum/Plasma	5.15

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	K151BDB-1	K111BDB-1
Tissue Culture	5 plates	K151BDB-2	K111BDB-2
Tissue Culture	25 plates	K151BDB-4	K111BDB-4
Ultra-Sensitive	1 plate	K151BDC-1	K111BDC-1
Ultra-Sensitive	5 plates	K151BDC-2	K111BDC-2
Ultra-Sensitive	25 plates	K151BDC-4	K111BDC-4
Base	25 plates	K151BDA-4	K111BDA-4
Base	250 plates	K151BDA-5	K111BDA-5