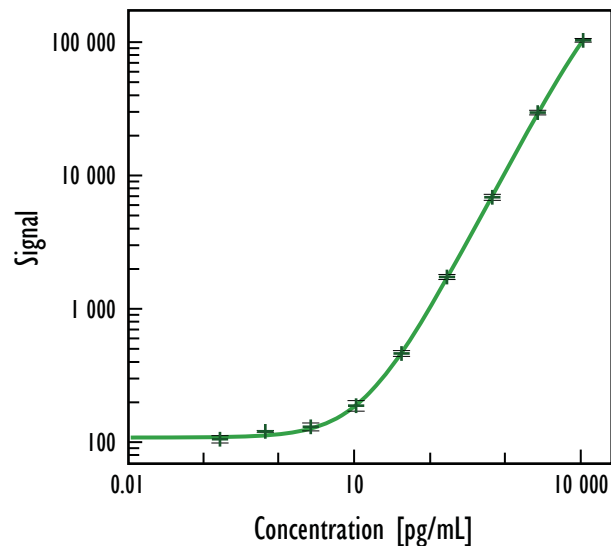


# MULTI-ARRAY<sup>®</sup> Human MIP-1 $\beta$ Tissue Culture Assay

## Detection of MIP-1 $\beta$ in Tissue Culture Samples

### Typical Standard Curve Data



Human MIP-1 $\beta$	
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	Catalog Number
Tissue Culture	1 plate	K151BDB-1
Tissue Culture	5 plates	K151BDB-2
Tissue Culture	25 plates	K151BDB-4
Ultra-Sensitive	1 plate	K151BDC-1
Ultra-Sensitive	5 plates	K151BDC-2
Ultra-Sensitive	25 plates	K151BDC-4

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

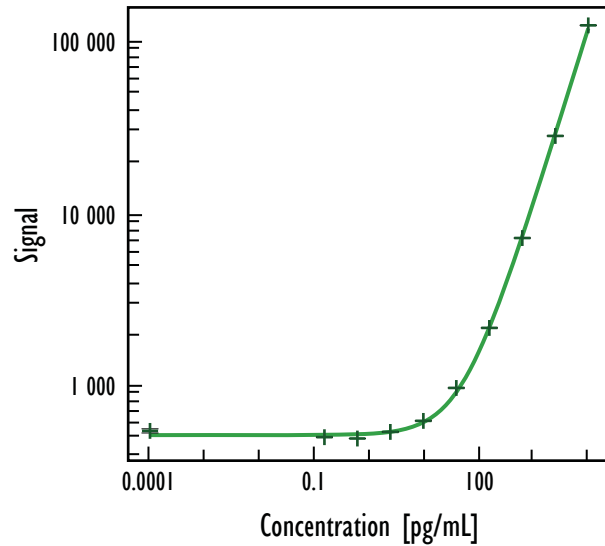


Meso Scale Discovery  
A division of Meso Scale Diagnostics, LLC.  
www.mesoscale.com

# MULTI-ARRAY<sup>®</sup> Human MIP-1 $\beta$ Ultra-Sensitive Assay

## Detection of MIP-1 $\beta$ in Human Serum and Plasma Samples

### Typical Standard Curve Data



Human MIP-1 $\beta$	
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	Catalog Number
Tissue Culture	1 plate	K151BDB-1
Tissue Culture	5 plates	K151BDB-2
Tissue Culture	25 plates	K151BDB-4
Ultra-Sensitive	1 plate	K151BDC-1
Ultra-Sensitive	5 plates	K151BDC-2
Ultra-Sensitive	25 plates	K151BDC-4