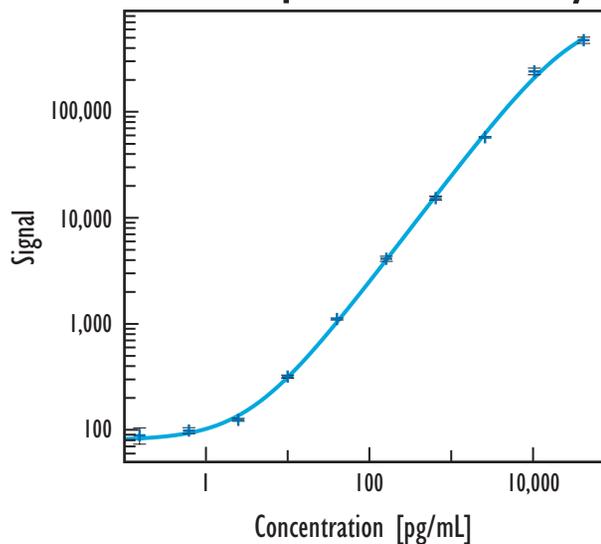


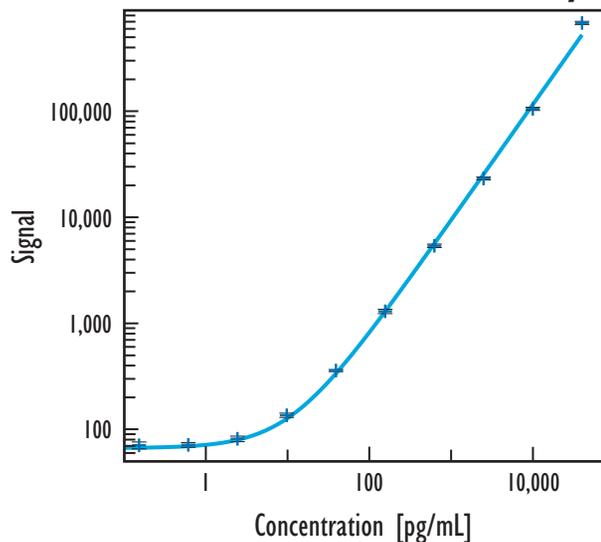
Mouse GM-CSF Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse GM-CSF		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	91	14.9
0.010	76	7.3
0.038	75	6.6
0.15	89	17.5
0.61	99	6.4
2.4	125	3.1
9.8	320	3.5
39	1,117	2.1
156	4,134	5.6
625	15,437	4.3
2,500	57,895	1.3
10,000	242,065	7.3
40,000	476,910	6.8

Serum / Plasma Assay



Mouse GM-CSF		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	59	9.8
0.15	71	7.5
0.61	72	5.2
2.4	82	6.1
9.8	136	4.7
39	359	2.3
156	1,298	4.8
625	5,390	3.8
2,500	23,243	3.0
10,000	105,475	3.5
40,000	680,775	3.4

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.8
Serum / Plasma	4.4

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	101
Serum	118
EDTA Plasma	113
Heparin Plasma	132

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=9)	9	n/d
EDTA Plasma (n=2)	1	n/d - 5.3
Heparin Plasma (n=9)	9	n/d

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.