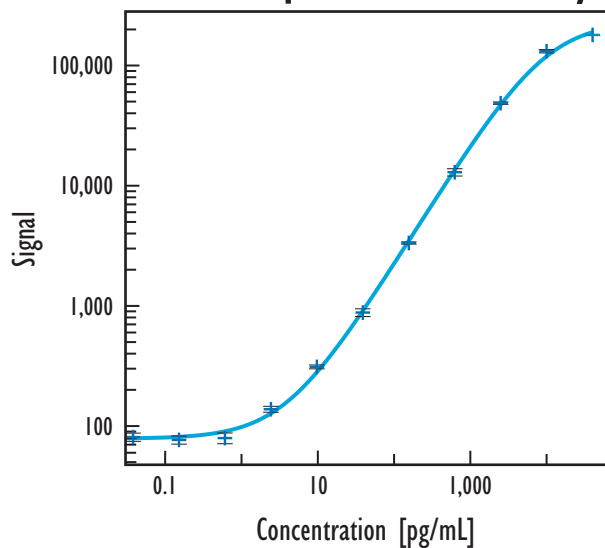


Mouse IL-1 β Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse IL-1 β		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	88	4.8
0.010	80	8.8
0.038	81	7.9
0.15	77	7.7
0.61	79	10.2
2.4	138	5.7
9.8	310	3.7
39	878	7.4
156	3,331	2.1
625	12,919	6.9
2,500	48,545	2.3
10,000	131,381	3.1
40,000	179,641	1.1

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.6
Serum / Plasma	1.5

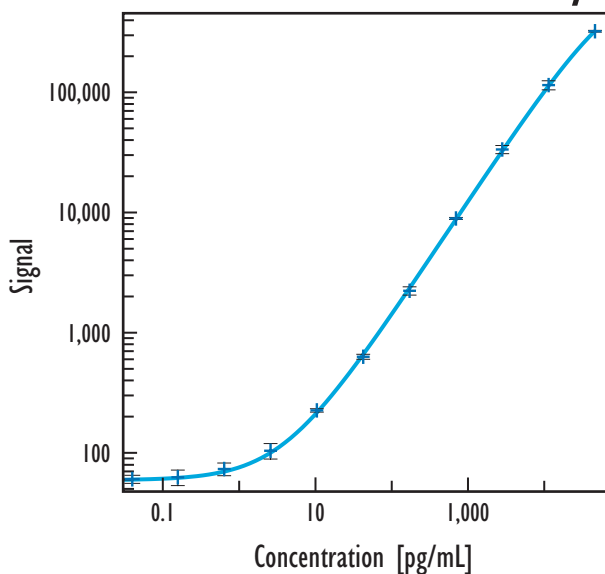
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	99
Serum	111
EDTA Plasma	101
Heparin Plasma	107

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.

Serum / Plasma Assay



Mouse IL-1 β		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	54	9.0
0.15	61	14.8
0.61	72	12.1
2.4	102	14.8
9.8	220	3.2
39	617	4.7
156	2,188	8.0
625	8,770	1.9
2,500	32,989	7.8
10,000	113,450	8.7
40,000	319,142	1.6

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	6	n/d - 2.1
Heparin Plasma (n=8)	6	n/d - 2.2

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.