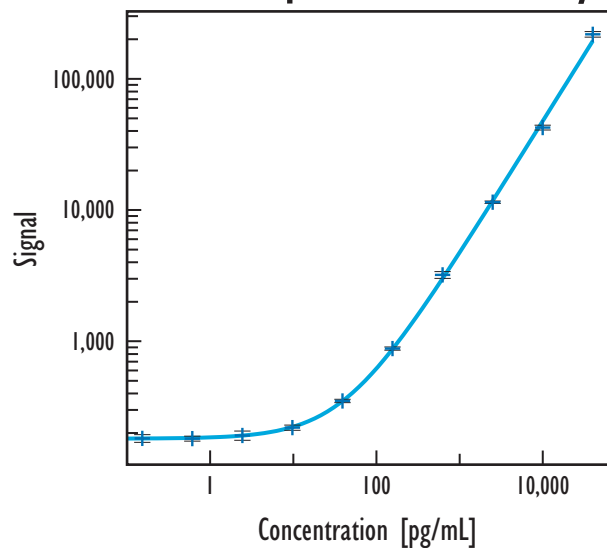


# Mouse IL-4 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

## Cell Supernatant Assay



Mouse IL-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	188	5.0
0.010	182	6.4
0.038	177	6.8
0.15	182	6.8
0.61	181	4.4
2.4	192	8.1
9.8	220	4.5
39	351	2.9
156	880	2.8
625	3,204	6.0
2,500	11,475	2.1
10,000	42,514	4.2
40,000	218,310	4.9

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	1.8
Serum / Plasma	3.5

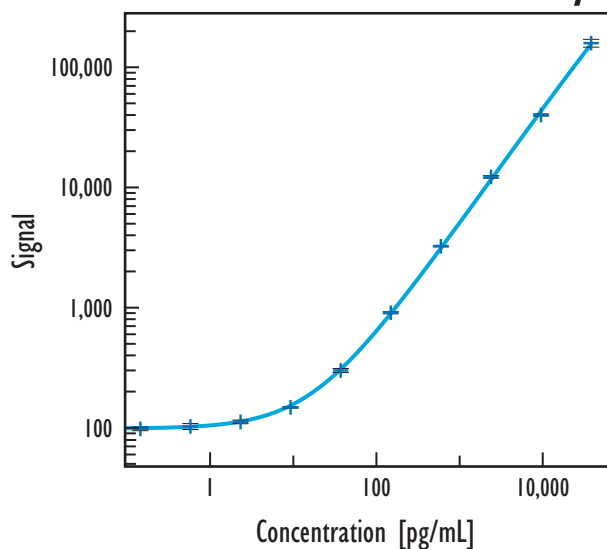
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

## Recoveries

Sample Type	Average % Recovery
Cell Supernatant	97
Serum	51
EDTA Plasma	54
Heparin Plasma	56

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-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	102	7.4
0.15	97	3.5
0.61	101	5.7
2.4	111	3.5
9.8	146	1.4
39	295	3.8
156	890	2.1
625	3,169	0.7
2,500	11,937	2.4
10,000	39,007	2.1
40,000	154,097	7.5

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	8	n/d
EDTA Plasma (n=1)	1	n/d
Heparin Plasma (n=8)	8	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.