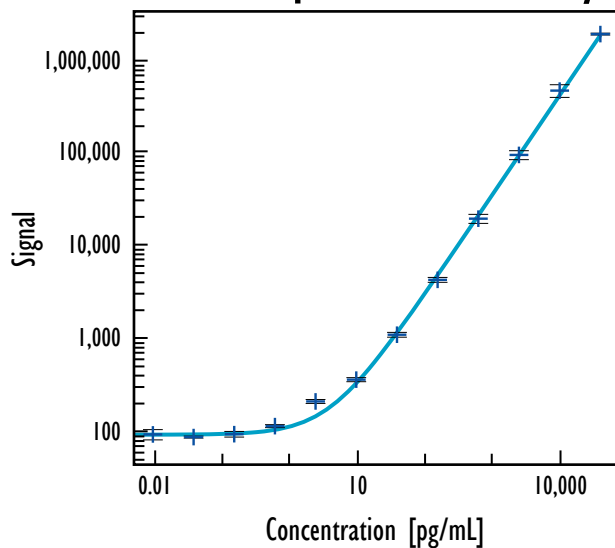


# Mouse IL-5 Cytokine Assay (MULTI-ARRAY® 96-Well Small Spot Plate)

## Cell Supernatant Assay



Mouse IL-5		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	80	8.2
0.010	91	12.7
0.038	85	2.4
0.15	91	7.0
0.61	111	3.7
2.4	206	4.9
9.8	355	5.1
39	1,085	5.9
156	4,257	6.1
625	19,528	11.2
2,500	95,506	11.1
10,000	474,267	15.7
40,000	1,937,253	2.3

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	1.0
Serum / Plasma	1.4

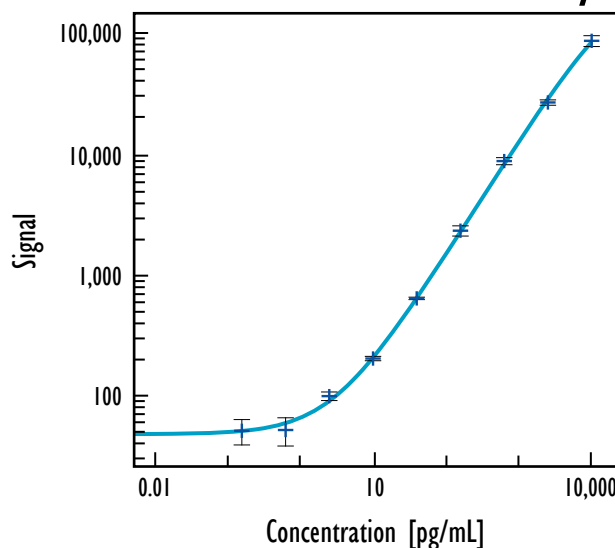
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

## Recoveries

Sample Type	Average % Recovery
Cell Supernatant	100
Serum	109
EDTA Plasma	119
Heparin Plasma	123

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-5		
Concentration (pg/mL)	Signal	
	Mean	%CV
0.001	54	22.9
0.15	52	23.7
0.61	53	26.2
2.4	101	8.1
9.8	205	4.1
39	644	2.3
156	2,331	9.7
625	8,763	6.8
2,500	26,682	5.1
10,000	86,067	10.3

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	0	3.1 - 15.2
Heparin Plasma (n=8)	1	1.7 - 12.8

Endogenous cytokine levels determined for different sample types across multiple samples.

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested.