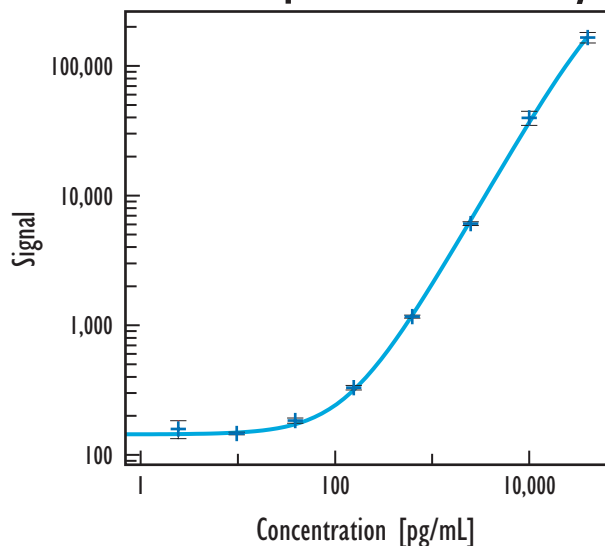


Mouse IL-10 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse IL-10		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	602	3.2
0.010	665	1.7
0.038	633	3.8
0.15	633	2.2
0.61	608	1.3
2.4	690	7.5
9.8	629	1.1
39	685	0.7
156	881	0.8
625	2,076	2.5
2,500	8,473	2.8
10,000	50,132	8.4
40,000	227,463	7.7

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	35
Serum / Plasma	42

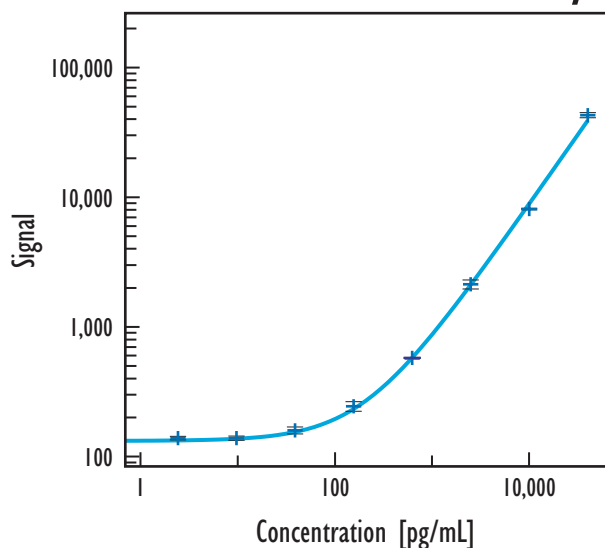
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	107
Serum	88
EDTA Plasma	97
Heparin Plasma	97

Spike recoveries were determined in each matrix over a range of spike levels. Each spike was tested in triplicate.

Serum / Plasma Assay



Mouse IL-10		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	132	5.9
0.15	128	6.2
0.61	131	4.7
2.4	141	3.1
9.8	141	3.7
39	162	6.2
156	248	8.5
625	584	1.0
2,500	2,156	7.9
10,000	8,186	1.3
40,000	43,330	4.5

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	8	n/d
Heparin Plasma (n=8)	7	n/d - 64

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.