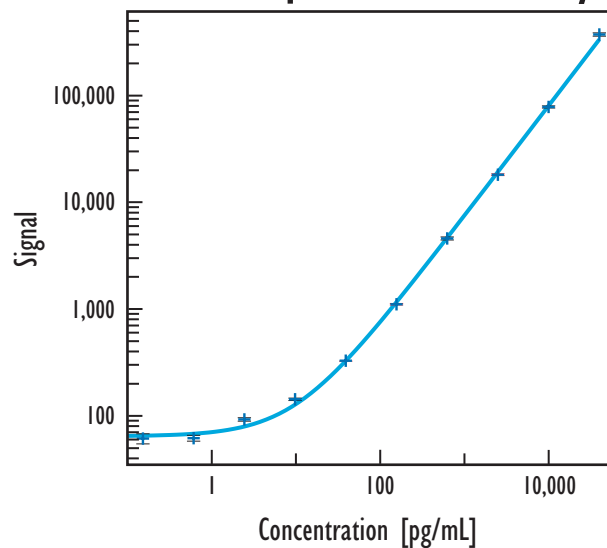


Mouse IL-2 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse IL-2		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	64	11.4
0.010	71	12.3
0.038	62	11.8
0.15	61	10.9
0.61	62	6.3
2.4	93	3.7
9.8	143	2.6
39	328	0.9
156	1,109	1.7
625	4,586	3.2
2,500	18,138	1.9
10,000	78,220	2.3
40,000	373,304	3.5

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	4.0
Serum / Plasma	10.4

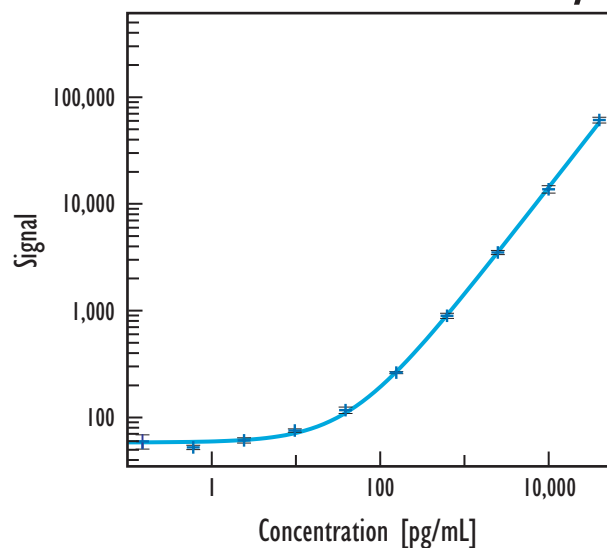
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	96
Serum	99
EDTA Plasma	88
Heparin Plasma	114

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-2		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	68	10.4
0.15	65	15.0
0.61	57	4.6
2.4	67	5.6
9.8	82	3.6
39	127	6.9
156	282	2.1
625	945	5.5
2,500	3,642	4.5
10,000	14,000	7.9
40,000	61,199	6.1

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	8	n/d
EDTA Plasma (n=2)	2	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.