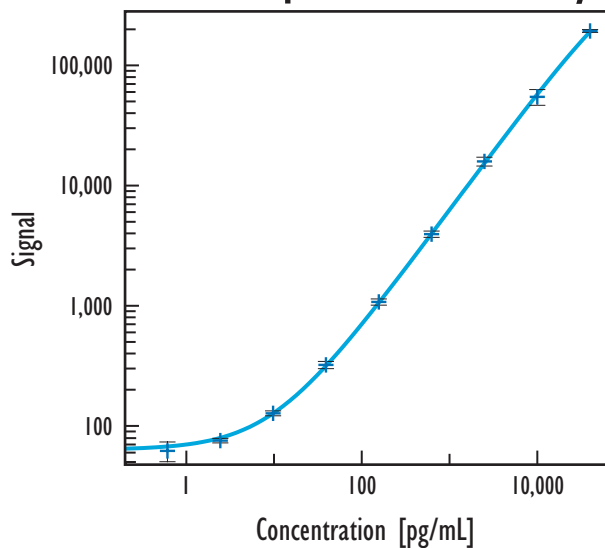


Mouse IL-12 Total Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse IL-12 (Total)		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	64	9.6
0.010	72	10.9
0.038	67	6.2
0.15	61	9.6
0.61	62	18.7
2.4	75	4.4
9.8	127	4.7
39	321	6.9
156	1,075	6.0
625	3,950	6.0
2,500	15,926	8.4
10,000	54,962	14.9
40,000	194,570	2.8

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	3.7
Serum / Plasma	6.9

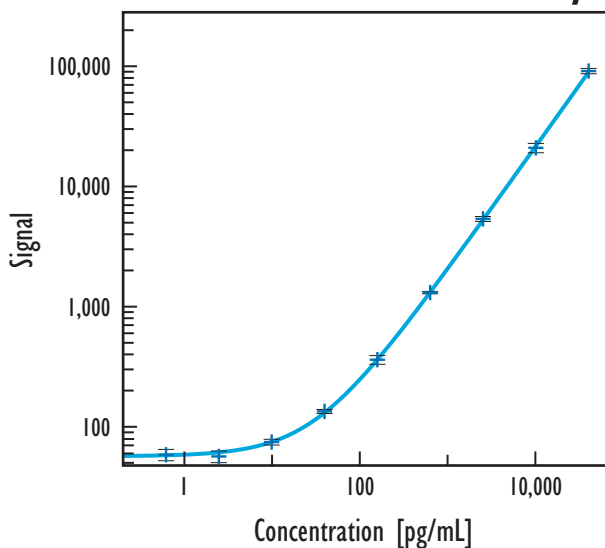
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	89
Serum	92
EDTA Plasma	97
Heparin Plasma	102

Spike recoveries were determined in cell culture over a range of spike levels from 625 to 10,000 pg/mL.

Serum / Plasma Assay



Mouse IL-12 (Total)		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	57	11.1
0.15	68	13.6
0.61	61	10.8
2.4	59	11.1
9.8	78	5.4
39	140	3.8
156	375	8.3
625	1,350	2.3
2,500	5,499	4.6
10,000	21,319	8.9
40,000	92,322	5.0

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	0	1,754 - 4,416
Heparin Plasma (n=8)	0	3,506 - 6,640

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.