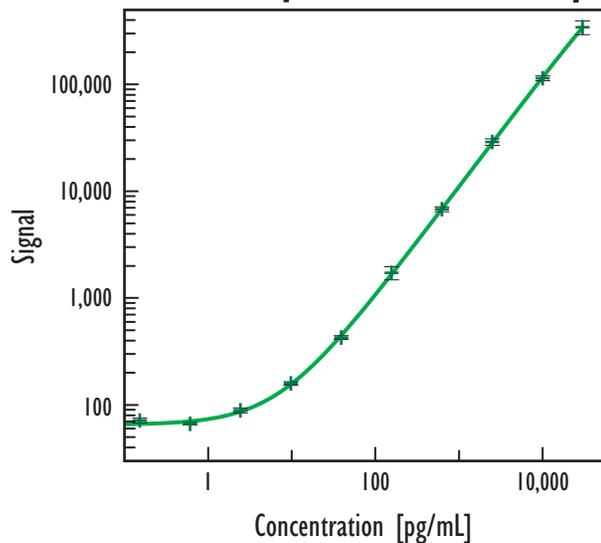


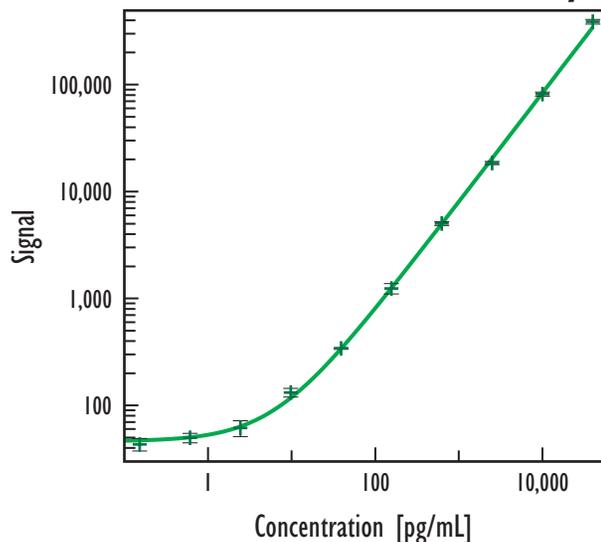
Human IL-12p70 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human IL-12p70		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	70	9.4
0.010	65	6.8
0.038	60	8.3
0.15	72	5.2
0.61	67	2.6
2.4	89	5.3
9.8	159	3.7
39	428	4.0
156	1,724	14.1
625	6,796	5.3
2,500	28,950	7.0
10,000	114,263	5.2
40,000	343,035	14.7

Serum / Plasma Assay



Human IL-12p70		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	51	14.0
0.15	43	13.4
0.61	50	10.2
2.4	62	16.8
9.8	132	9.3
39	341	1.6
156	1,240	11.1
625	5,027	4.1
2,500	18,497	3.4
10,000	81,654	4.4
40,000	387,908	5.0

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	2.8
Serum	3.5
Plasma	8.8

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	99
Serum	110
EDTA Plasma	95
Heparin Plasma	100

Spike recoveries were determined in each matrix over a range of spike levels from 9.8 to 313 pg/mL. Each spike was tested in ≥ 3 replicates.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=5)	5	n/d
Heparin Plasma (n=5)	5	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.