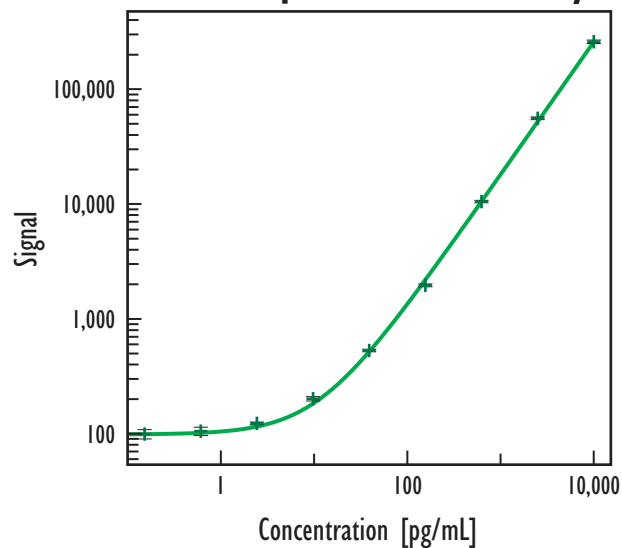


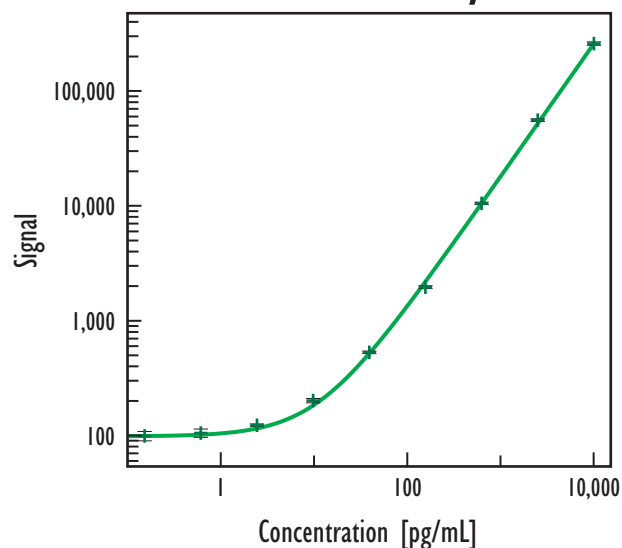
Human IL-4 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human IL-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	60	13.3
0.010	60	8.5
0.038	54	10.8
0.15	50	15.0
0.61	49	14.5
2.4	77	9.2
9.8	105	7.4
39	258	4.6
156	878	1.8
625	3,811	6.6
2,500	15,077	8.0
10,000	59,745	7.8
40,000	323,509	2.2

Serum Assay



Human IL-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	55	10.0
0.15	49	8.3
0.61	52	14.0
2.4	78	8.6
9.8	129	4.9
39	327	3.3
156	1,055	3.9
625	3,861	7.0
2,500	14,823	8.9
10,000	58,352	2.3
40,000	271,778	10.3

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	5.3
Serum	3.6
Plasma	12.0

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	102
Serum	81
EDTA Plasma	111
Heparin Plasma	133

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.