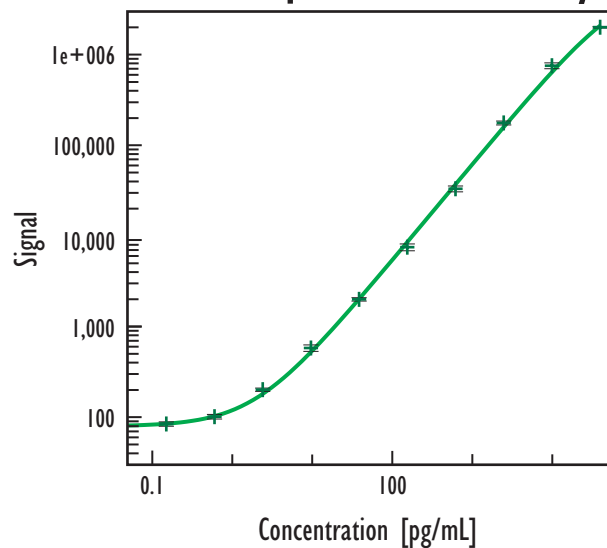


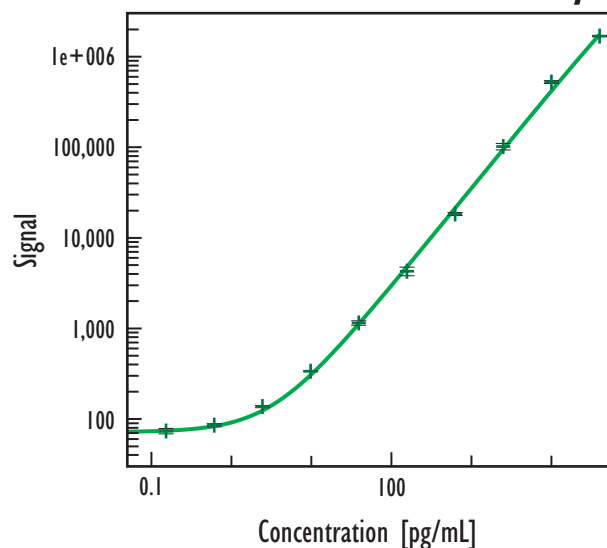
Human IL-5 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human IL-5		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	78	9.9
0.010	79	2.2
0.038	76	5.1
0.15	83	5.5
0.61	100	5.9
2.4	198	4.2
9.8	571	8.2
39	1,965	4.3
156	7,378	8.6
625	32,492	7.4
2,500	172,806	4.9
10,000	739,044	7.4
40,000	1,961,628	2.0

Serum / Plasma Assay



Human IL-5		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	67	8.1
0.15	74	6.9
0.61	85	4.0
2.4	137	2.7
9.8	336	1.2
39	1,142	5.5
156	4,252	10.8
625	18,181	3.8
2,500	100,642	8.3
10,000	517,480	3.6
40,000	1,663,160	1.5

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.7
Serum	1.2
Plasma	3.4

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	100
Serum	91
EDTA Plasma	93
Heparin Plasma	106

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=4)	4	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.