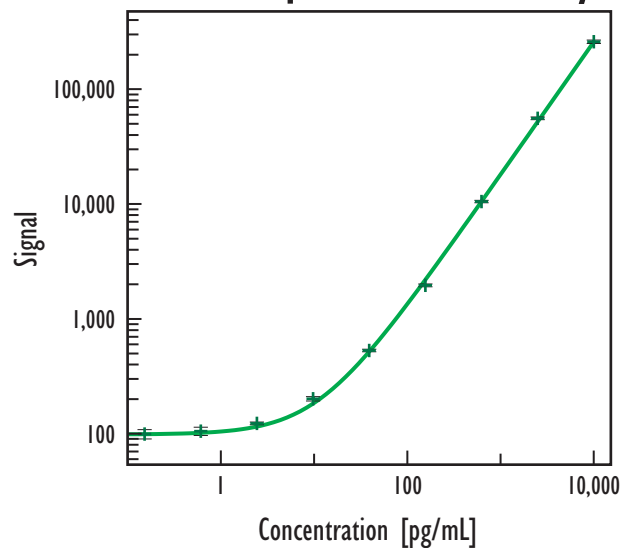


Human IL-10 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



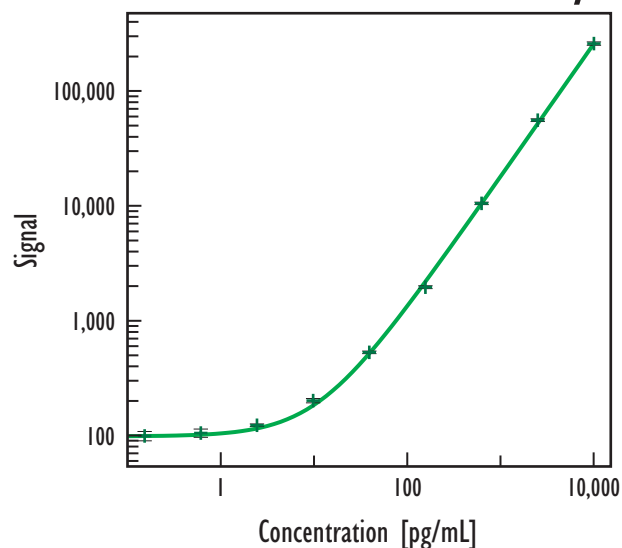
Human IL-10		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	72	8.1
0.010	60	6.2
0.038	66	9.3
0.15	63	12.1
0.61	60	7.8
2.4	73	6.2
9.8	101	5.2
39	187	3.4
156	590	1.8
625	2,067	11.8
2,500	10,250	3.7
10,000	44,149	5.2
40,000	188,697	8.8

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	9.1
Serum	23.4
Plasma	36.6

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

Serum / Plasma Assay



Human IL-10		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	62	9.3
0.15	56	8.1
0.61	55	10.0
2.4	56	6.7
9.8	79	5.2
39	109	7.5
156	252	2.5
625	842	3.8
2,500	3,248	10.0
10,000	13,317	8.4
40,000	77,477	9.1

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	9	n/d
EDTA Plasma (n=5)	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.