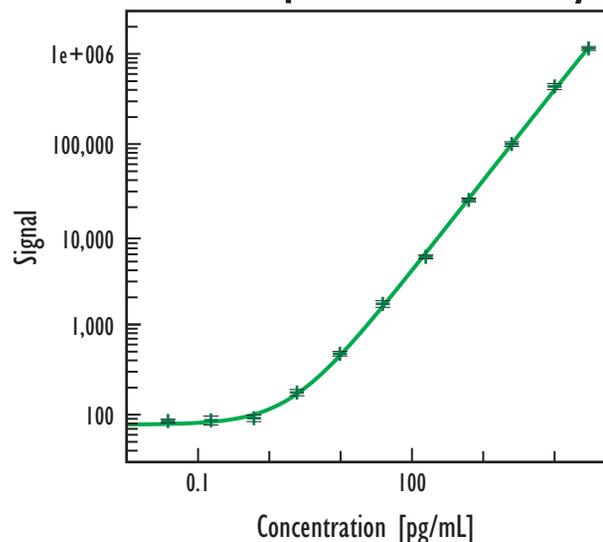


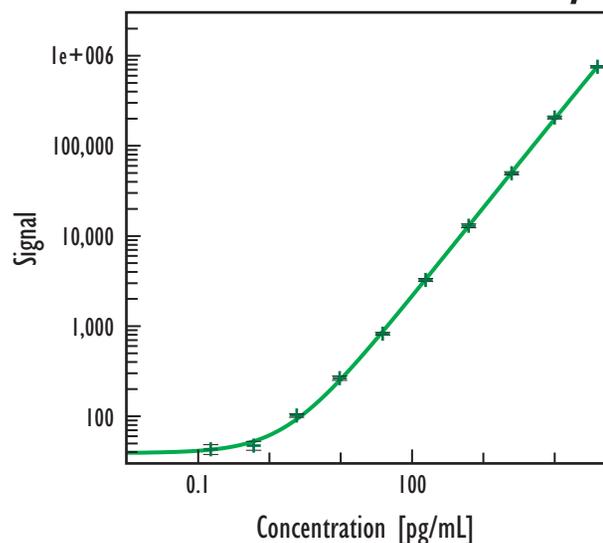
Human TNF- α Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human TNF- α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	75	9.1
0.010	77	5.5
0.038	85	5.0
0.15	87	11.3
0.61	92	8.3
2.4	176	8.0
9.8	475	5.9
39	1,691	8.3
156	5,615	4.8
625	23,992	4.7
2,500	99,871	6.0
10,000	433,810	7.6
40,000	1,144,978	5.0

Serum / Plasma Assay



Human TNF- α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	41	17.5
0.15	43	12.3
0.61	47	11.2
2.4	102	4.5
9.8	264	5.6
39	821	3.4
156	3,235	3.9
625	12,891	4.9
2,500	49,351	4.0
10,000	204,799	3.8
40,000	746,301	3.0

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.7
Serum	1.1
Plasma	2.7

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	106
Serum	90
EDTA Plasma	102
Heparin Plasma	111

Spike recoveries were determined in each matrix over a range of spike levels from 9.8 to 625 pg/mL. Each spike was tested in ≥ 3 replicates.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	0	3.3 - 9.3
EDTA Plasma (n=5)	4	n/d - 2.9
Heparin Plasma (n=5)	1	n/d - 4.1

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.