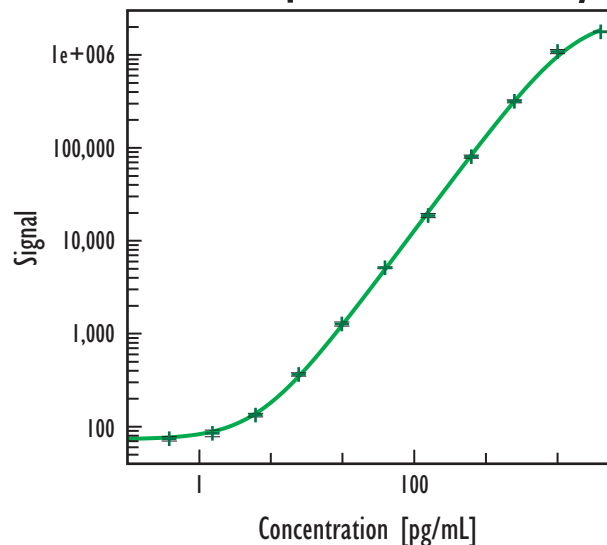


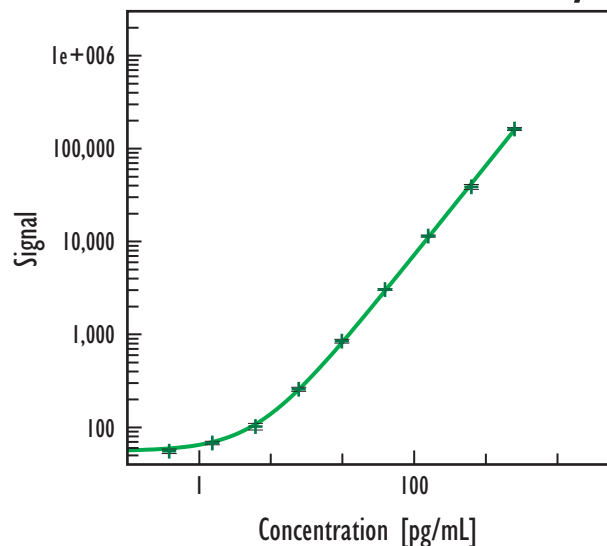
Human TNF- α Ultrasensitive Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human TNF- α US		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	81	7.4
0.010	74	4.5
0.038	74	5.5
0.15	85	8.2
0.61	134	4.4
2.4	365	4.2
9.8	1,275	5.1
39	5,130	2.1
156	18,732	5.0
625	80,597	3.9
2,500	317,523	3.7
10,000	1,089,784	5.1
40,000	1,776,323	1.6

Serum / Plasma Assay



Human TNF- α US		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	61	10.0
0.15	59	11.3
0.61	56	5.6
2.4	68	4.3
9.8	102	8.2
39	258	5.0
156	847	4.7
625	3,036	2.1
2,500	11,396	2.5
10,000	38,595	6.0
40,000	161,589	3.5

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.2
Serum	0.3
Plasma	0.7

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	100
Serum	88
EDTA Plasma	106
Heparin Plasma	114

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)	0	1.9 - 5.3
EDTA Plasma (n=5)	1	n/d - 2.2
Heparin Plasma (n=5)	0	1.6 - 3.8

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.