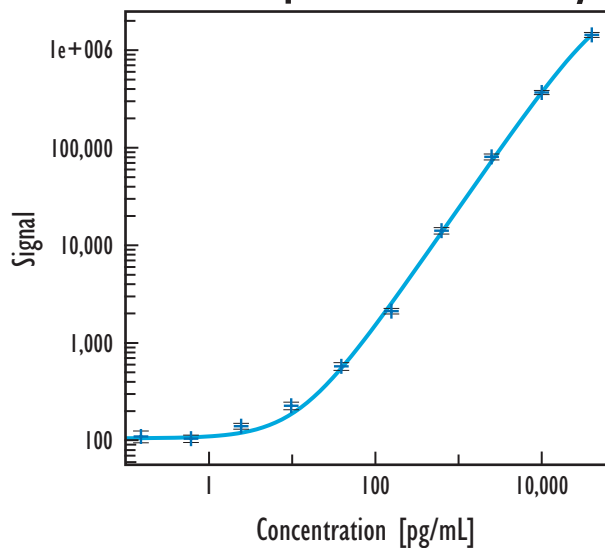


Mouse TNF- α Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse TNF- α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	99	6.3
0.010	101	7.4
0.038	99	5.2
0.15	110	13.8
0.61	105	8.6
2.4	140	6.9
9.8	227	8.9
39	576	9.0
156	2,120	6.5
625	14,129	7.5
2,500	80,739	7.0
10,000	368,869	4.9
40,000	1,434,346	5.8

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	3.7
Serum / Plasma	3.4

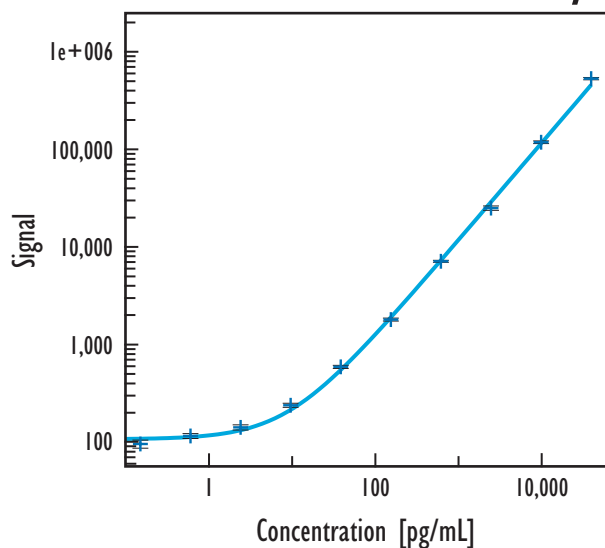
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	95
Serum	117
EDTA Plasma	121
Heparin Plasma	103

Spike recoveries were determined in each matrix over a range of spike levels from 39 to 625 pg/mL. Each spike was tested in triplicate.

Serum / Plasma Assay



Mouse TNF- α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	72	6.9
0.15	62	9.9
0.61	76	6.0
2.4	94	6.4
9.8	162	5.0
39	418	3.9
156	1,346	3.7
625	5,681	2.4
2,500	21,174	5.3
10,000	107,843	3.2
40,000	519,366	2.9

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=9)	8	n/d - 5.6
EDTA Plasma (n=2)	1	n/d - 12
Heparin Plasma (n=9)	9	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.