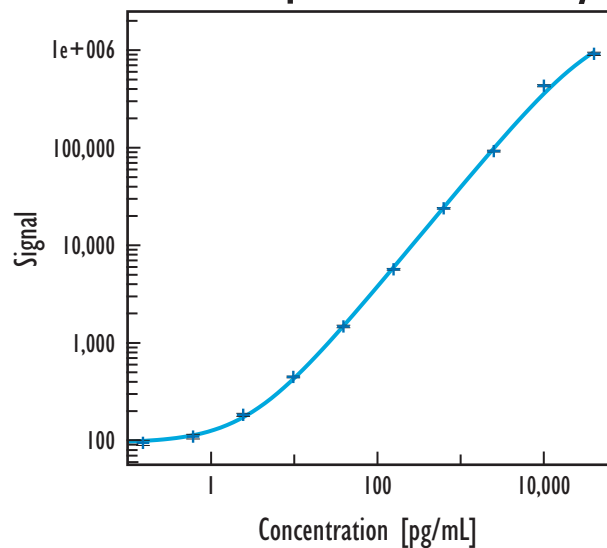


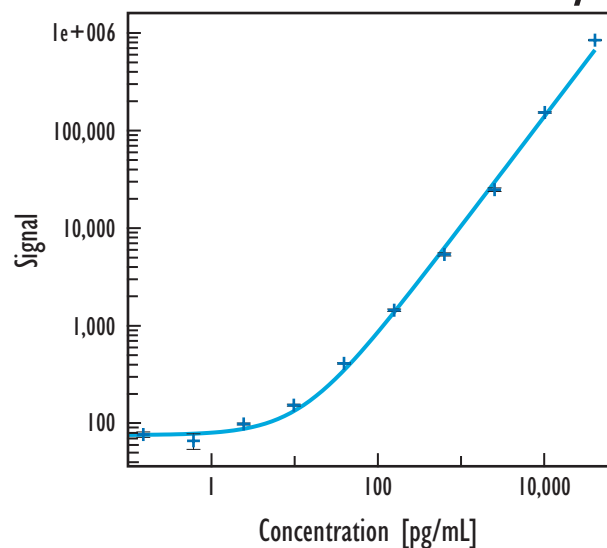
Mouse IFN- γ Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse IFN- γ		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	87	7.5
0.010	104	3.6
0.038	95	8.9
0.15	93	6.2
0.61	108	5.6
2.4	181	3.7
9.8	447	2.0
39	1,493	3.0
156	5,841	2.0
625	25,073	1.9
2,500	98,187	2.0
10,000	467,557	2.5
40,000	1,001,449	3.9

Serum / Plasma Assay



Mouse IFN- γ		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	77	11.1
0.15	76	6.5
0.61	65	18.2
2.4	97	2.3
9.8	152	2.3
39	407	1.6
156	1,431	2.8
625	5,352	3.7
2,500	24,722	4.2
10,000	153,251	1.7
40,000	847,004	1.0

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.8
Serum / Plasma	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	114
EDTA Plasma	113
Heparin Plasma	121

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.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	0	27 - 70
Heparin Plasma (n=8)	0	12 - 60

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.