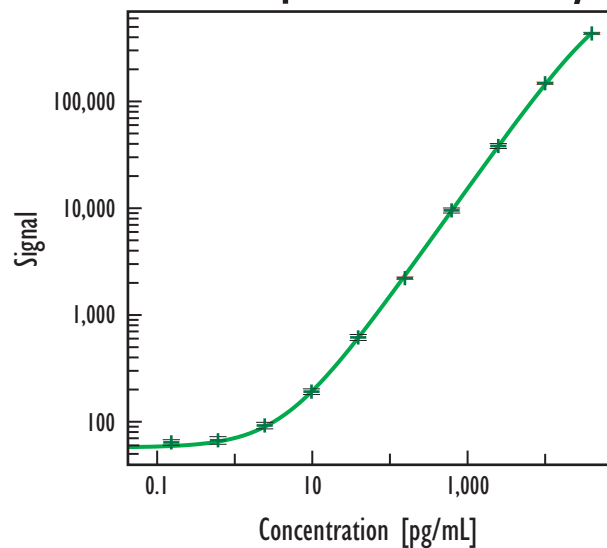


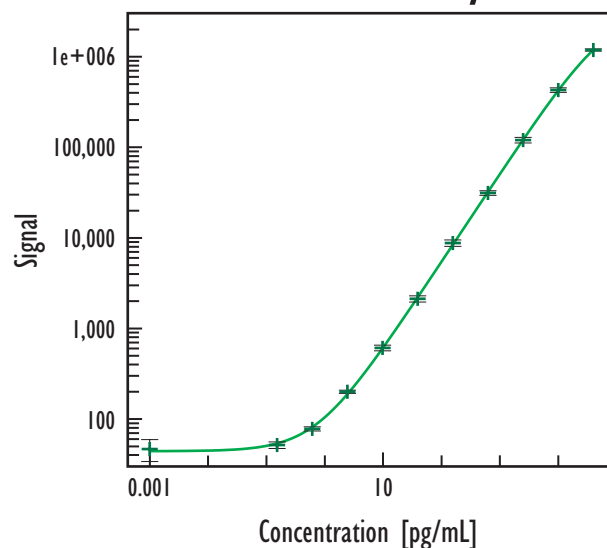
# Human IFN- $\gamma$ Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

## Cell Supernatant Assay



Human IFN- $\gamma$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	52	7.3
0.010	56	13.9
0.038	60	5.8
0.15	64	6.0
0.61	67	8.2
2.4	92	7.0
9.8	192	6.1
39	617	6.5
156	2,213	2.6
625	9,531	5.5
2,500	38,273	5.3
10,000	148,033	2.5
40,000	433,411	2.5

## Serum Assay



Human IFN- $\gamma$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	54	7.7
0.15	66	7.5
0.61	83	1.4
2.4	155	12.0
9.8	412	7.8
39	1,300	3.4
156	5,113	5.0
625	19,821	4.9
2,500	84,323	2.4
10,000	284,282	11.3
40,000	662,964	2.7

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	1.5
Serum	4.4
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	91
Serum	92
EDTA Plasma	106
Heparin Plasma	107

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  $\geq 3$  replicates.

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	8	n/d
EDTA Plasma (n=4)	4	n/d
Heparin Plasma (n=4)	3	n/d - 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.