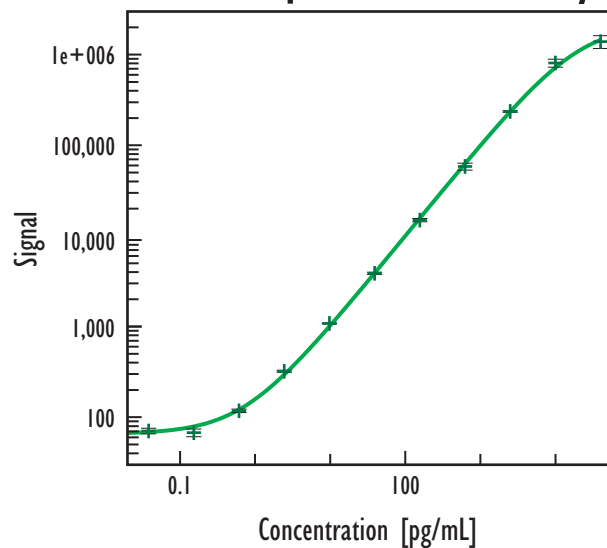


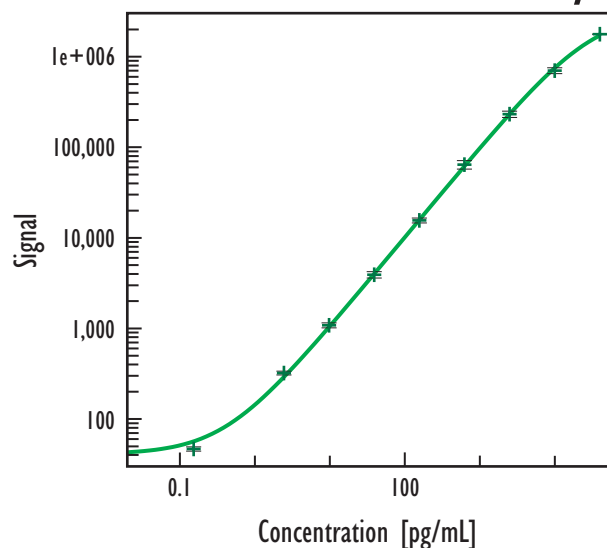
Human IL-8 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human IL-8		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	78	14.7
0.010	67	11.7
0.038	71	7.1
0.15	68	9.7
0.61	118	4.4
2.4	322	2.5
9.8	1,083	1.1
39	3,869	2.9
156	14,962	4.2
625	58,406	8.8
2,500	236,433	2.2
10,000	809,925	10.1
40,000	1,394,565	16.1

Serum / Plasma Assay



Human IL-8		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	53	9.4
0.15	47	5.3
2.4	322	4.9
9.8	1,083	6.3
39	3,915	8.1
156	15,579	6.2
625	64,351	10.7
2,500	231,900	8.0
10,000	701,686	7.4
40,000	1,771,058	0.6

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.3
Serum	0.3
Plasma	0.6

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	90
Serum	87
EDTA Plasma	85
Heparin Plasma	93

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=9)	0	4.2 - 42.6
EDTA Plasma (n=5)	0	0.9 - 2.1
Heparin Plasma (n=5)	0	1.8 - 5.2

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.