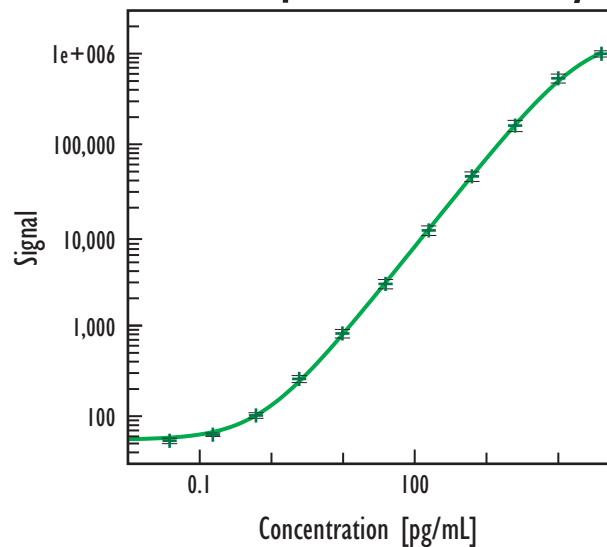


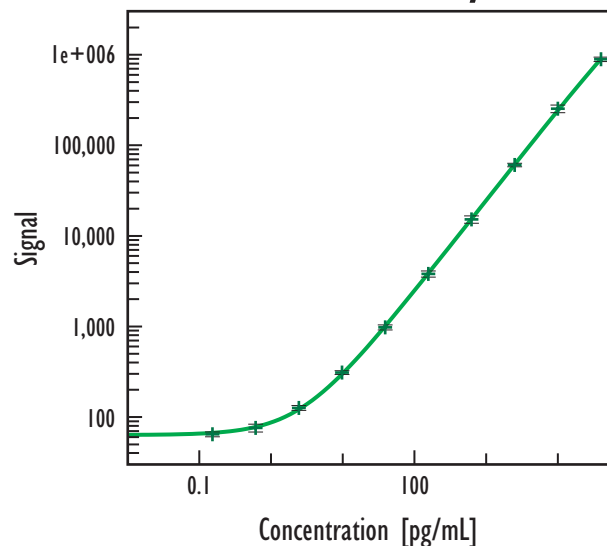
Human IL-1 β Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Human IL-1 β		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	65	8.0
0.010	56	20.3
0.038	54	7.0
0.15	63	4.2
0.61	102	6.9
2.4	258	8.8
9.8	819	10.8
39	2,877	11.8
156	11,188	12.2
625	44,110	12.0
2,500	159,937	14.2
10,000	531,024	11.3
40,000	988,914	7.8

Serum Assay



Human IL-1 β		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	65	7.9
0.15	64	6.1
0.61	75	10.0
2.4	125	6.1
9.8	308	4.7
39	971	6.4
156	3,781	7.8
625	15,147	9.4
2,500	60,426	4.4
10,000	252,960	9.8
40,000	889,148	5.6

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	0.3
Serum	2.4
Plasma	2.6

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	99
Serum	81
EDTA Plasma	89
Heparin Plasma	107

Spike recoveries were determined in each matrix over a range of spike levels from 9.8 to 625 pg/mL. Each spike was tested in ≥ 3 replicates.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=4)	4	n/d
Heparin Plasma (n=5)	5	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.