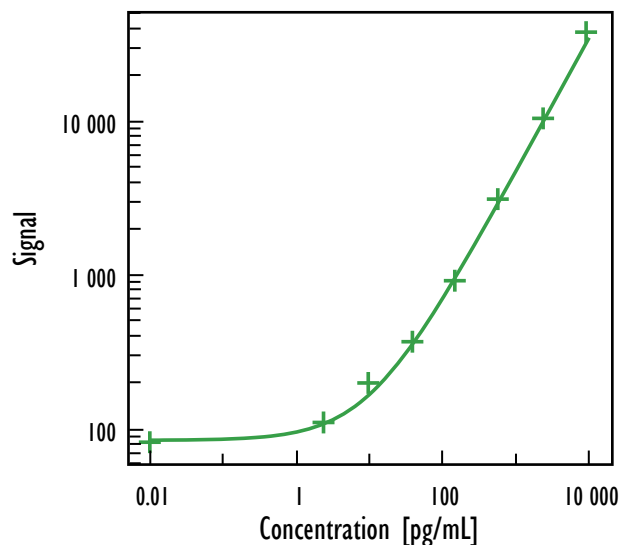


# MULTI-ARRAY<sup>®</sup> Human IL-12 Total Cytokine Assay

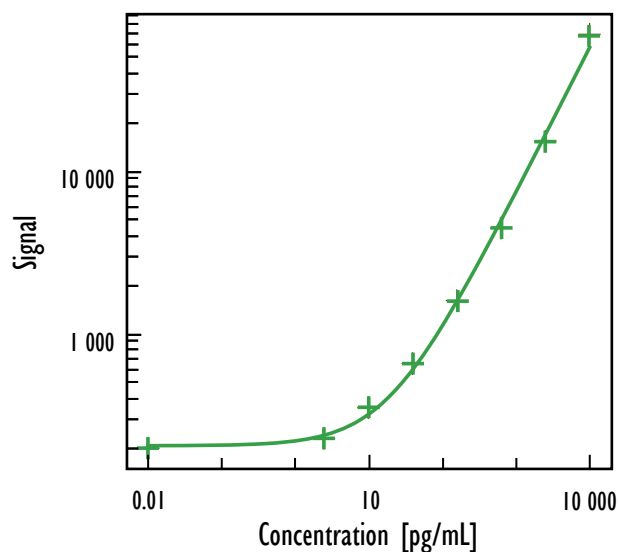
## Tissue Culture Kit



Human IL-12 Total		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	80	3.8
2.4	110	5.5
9.8	195	2.7
39	365	7.9
156	910	3.9
625	3020	11.5
2500	10089	7.5
10000	37443	9.8

Protocol	Detection Limit (pg/mL)
Tissue Culture	2.5

## Ultra-Sensitive Kit



Human IL-12 Total		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	198	3.0
2.4	228	4.3
9.8	352	3.1
39	657	7.1
156	1608	6.9
625	4500	10.3
2500	15429	1.3
10000	70482	3.8

Protocol	Detection Limit (pg/mL)
Serum/Plasma	2.0

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

## Endogenous Levels

- Pooled normal human serum and heparin plasma samples were measured.

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=2)	0	38 - 49
Heparin Plasma (n=2)	0	38 - 157

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K15IAPB-1
Tissue Culture	5 plates	K15IAPB-2
Tissue Culture	25 plates	K15IAPB-4
Ultra-Sensitive	1 plate	K15IAPC-1
Ultra-Sensitive	5 plates	K15IAPC-2
Ultra-Sensitive	25 plates	K15IAPC-4
Base	25 plates	K15IAPA-4
Base	250 plates	K15IAPA-5

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested.

# MULTI-ARRAY<sup>®</sup> Human IL-12 Total Cytokine Assay

## Dilution Linearity

- Human serum and heparin plasma pooled samples were evaluated for linear dilution. Pooled samples were spiked with calibrator and then diluted with Diluent 2. Recovery is calculated from the measured concentration of the previous dilution.

$$\% \text{ Recovery} = (\text{measured} \times \text{dilution factor}) / \text{expected} \times 100$$

	Human IL-12 Total			
	Fold Dilution	Conc. (pg/mL)	% CV	% Recovery
Serum	1	2762	3.4	
	2	2878	7.3	104
	4	3140	3.9	109
	8	2848	0.6	91
Heparin Plasma	1	2472	4.7	
	2	2752	5.0	111
	4	3000	2.3	109
	8	3048	0.1	102

## Spike Recovery

- Human serum, heparin plasma, and EDTA plasma pooled samples were spiked with calibrators at multiple levels throughout the range of the assay. Each spike was tested in  $\geq 3$  replicates.

$$\% \text{ Recovery} = \text{measured} / \text{expected} \times 100$$

Human IL-12 Total	Spike Level (pg/mL)	Measured Conc. (pg/mL)	% CV	% Recovery
Serum	0	38	5.5	
	31.5	69	2.5	99
	295	316	4.2	95
	3207	2998	2.1	92
EDTA Plasma	0	8	1.4	
	31.5	39	3.5	99
	295	247	3.7	82
	3207	2682	1.0	83
Heparin Plasma	0	157	2.7	
	31.5	191	6.2	101
	295	452	0.6	100
	3207	3495	9.0	104