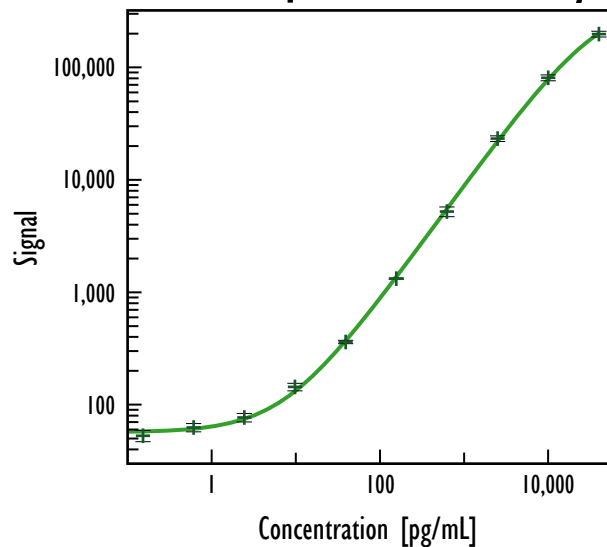


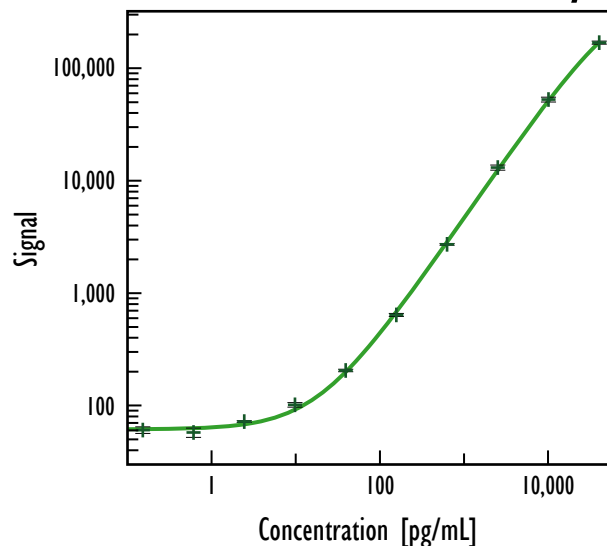
# Human IL-13 Cytokine Assay (MULTI-ARRAY<sup>®</sup> 96-Well 4 Spot Plate)

## Cell Supernatant Assay



Human IL-13		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	53	11.8
0.010	61	2.8
0.038	60	13.8
0.15	53	11.4
0.61	63	8.5
2.4	77	8.7
9.8	144	7.7
39	362	3.4
156	1,323	1.3
625	5,228	9.9
2,500	23,295	5.7
10,000	80,950	6.0
40,000	198,129	5.9

## Serum / Plasma Assay



Human IL-13		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	62	10.3
0.15	61	6.8
0.61	58	9.5
2.4	72	1.7
9.8	102	4.7
39	206	2.6
156	641	3
625	2,716	1.7
2,500	13,069	5.2
10,000	52,568	4.9
40,000	168,508	3.1

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	3.4
Serum	8.1
Plasma	18.3

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

## Recoveries

Sample Type	Average % Recovery
Cell Supernatant	96
Serum	79
EDTA Plasma	121
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=10)	10	n/d
EDTA Plasma (n=5)	5	n/d
Heparin Plasma (n=5)	4	n/d - 39.2

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.



Meso Scale Discovery<sup>®</sup>  
A division of Meso Scale Diagnostics, LLC.  
www.mesoscale.com<sup>™</sup>