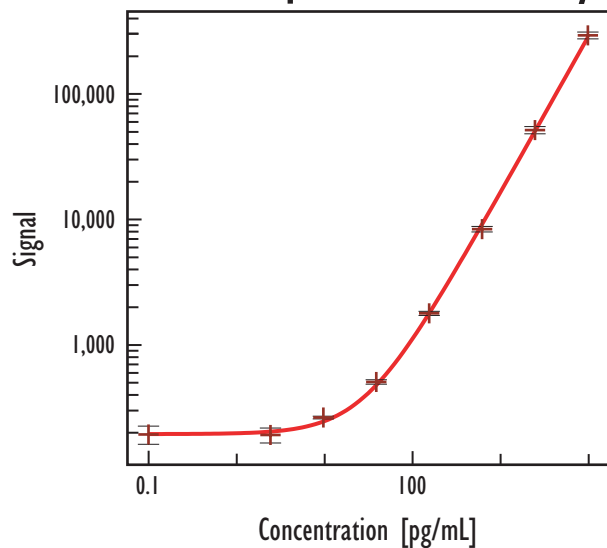


Rat IL-4 Cytokine Assay (MULTI-SPOT[®] 96-Well 7-Spot Plate)

Cell Supernatant Assay



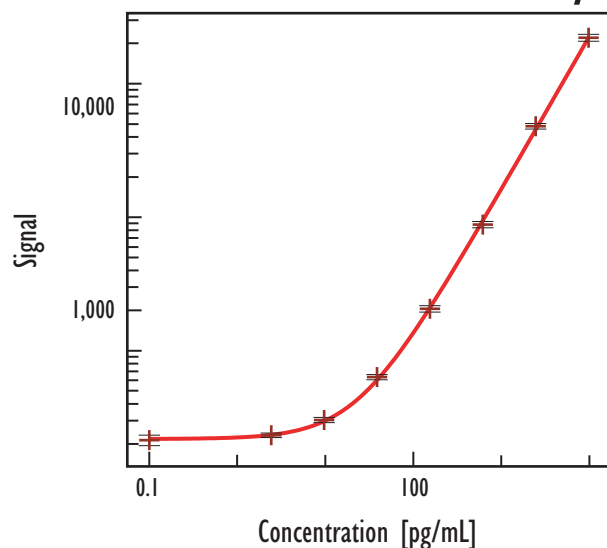
Rat IL-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	194	16.5
2.4	192	13.7
9.8	264	2.5
39	508	4.0
156	1,788	3.8
625	8,384	5.0
2,500	51,605	6.7
10,000	293,341	6.3

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	12.6
Serum / Plasma	7.5

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

Serum / Plasma Assay



Rat IL-4		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	214	9.1
2.4	233	3.9
9.8	303	4.5
39	635	4.5
156	2,061	5.5
625	8,816	5.4
2,500	48,109	4.7
10,000	221,624	6.0

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	0	10.2 - 21.8
Heparin Plasma (n=10)	4	n/d - 19.6

Endogenous cytokine levels (pg/mL) were determined for different sample types across multiple samples. For each sample type, samples from 5 male and 5 female rats were tested. Serum, EDTA plasma, and heparin plasma samples were not obtained from the same animal. (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.