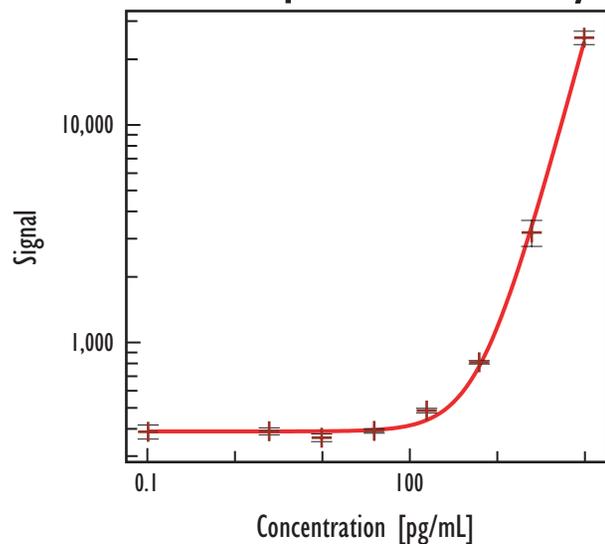


Rat MIP-3 α Cytokine Assay (MULTI-SPOT[®] 96-Well 7-Spot Plate)

Cell Supernatant Assay



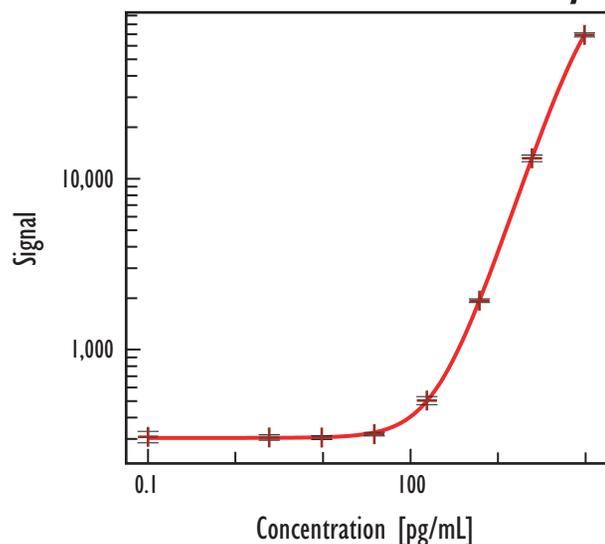
Rat MIP-3 α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	388	7.4
2.4	390	3.7
9.8	365	4.1
39	392	2.4
156	486	2.4
625	810	1.8
2,500	3,200	13.7
10,000	25,176	7.1

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	190.3
Serum / Plasma	57.9

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

Serum / Plasma Assay



Rat MIP-3 α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	309	7.7
2.4	306	3.7
9.8	306	2.7
39	320	2.4
156	504	5.5
625	1,935	2.5
2,500	13,170	4.4
10,000	69,385	2.7

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	8	n/d - 88.1
EDTA Plasma (n=10)	4	n/d - 425.3
Heparin Plasma (n=10)	6	n/d - 55.9

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.