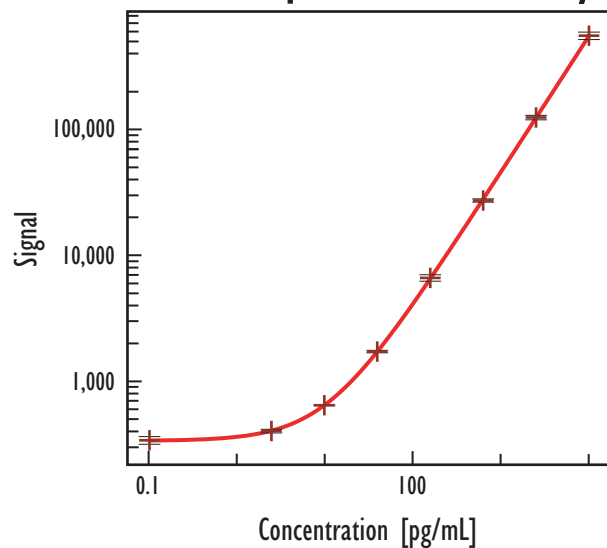


Rat GM-CSF Cytokine Assay (MULTI-SPOT® 96-Well 7-Spot Plate)

Cell Supernatant Assay



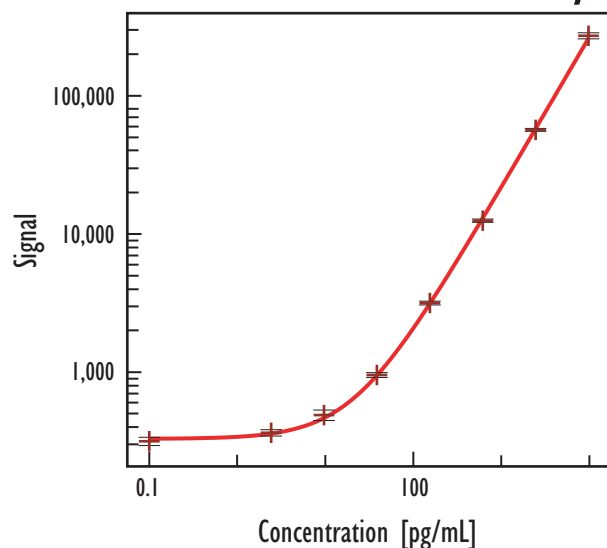
Rat GM-CSF		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	340	7.0
2.4	402	3.3
9.8	645	0.7
39	1,723	2.2
156	6,627	5.9
625	27,313	3.0
2,500	124,580	4.1
10,000	556,204	6.7

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	2.8
Serum / Plasma	4.7

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

Serum / Plasma Assay



Rat GM-CSF		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	316	6.8
2.4	364	5.2
9.8	490	8.7
39	955	4.0
156	3,170	3.2
625	12,514	2.7
2,500	56,884	2.3
10,000	272,992	5.0

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	4	n/d - 13.9
Heparin Plasma (n=10)	9	n/d - 7.4

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.