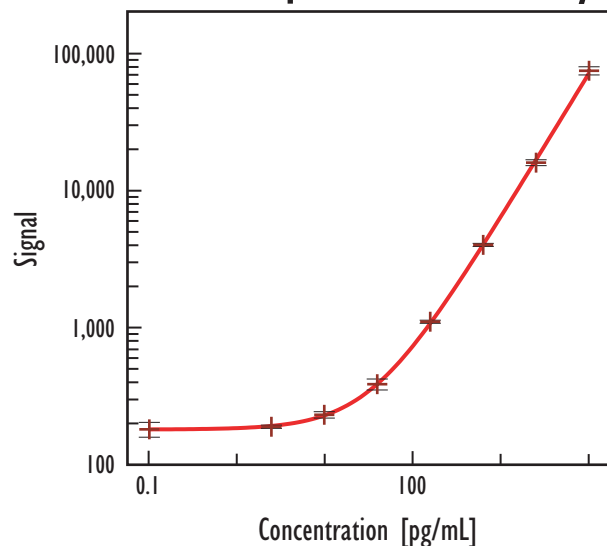


# Rat TNF- $\alpha$ Cytokine Assay (MULTI-SPOT<sup>®</sup> 96-Well 7-Spot Plate)

## Cell Supernatant Assay



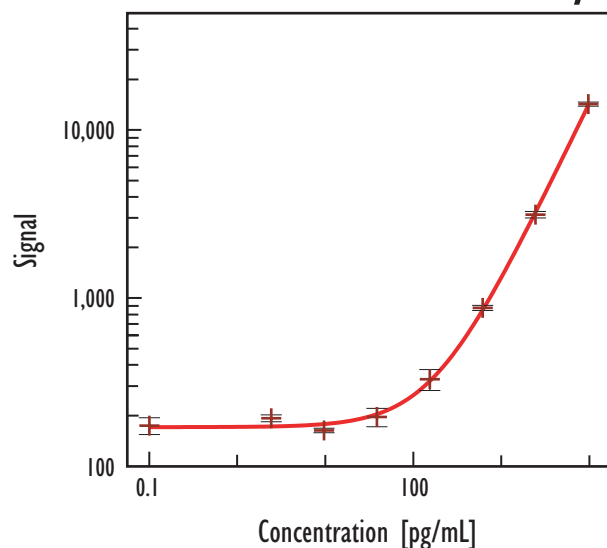
Rat TNF- $\alpha$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	181	12.4
2.4	189	2.6
9.8	231	5.3
39	386	9.1
156	1,103	2.6
625	4,002	2.6
2,500	15,927	5.0
10,000	74,420	6.9

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	18.9
Serum / Plasma	63.2

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

## Serum / Plasma Assay



Rat TNF- $\alpha$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	174	11.4
2.4	193	4.7
9.8	163	3.1
39	196	12.5
156	329	14.3
625	874	3.5
2,500	3,135	4.5
10,000	14,256	2.8

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	0	79.4 - 254.7
Heparin Plasma (n=10)	0	55.5 - 137.7

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.