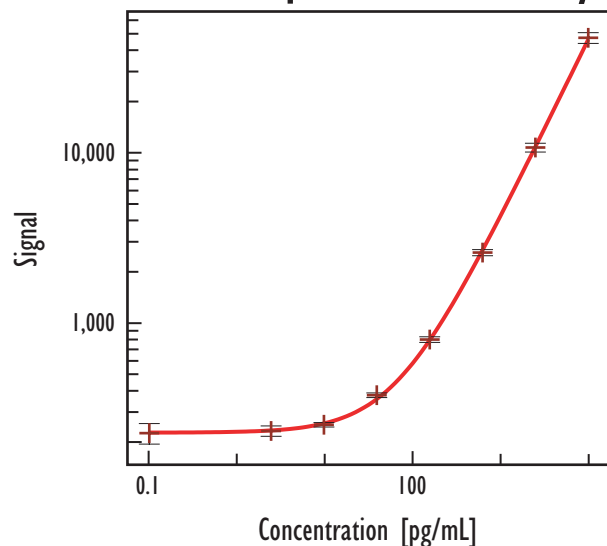


# Rat IL-1 $\beta$ Cytokine Assay (MULTI-SPOT<sup>®</sup> 96-Well 7-Spot Plate)

## Cell Supernatant Assay



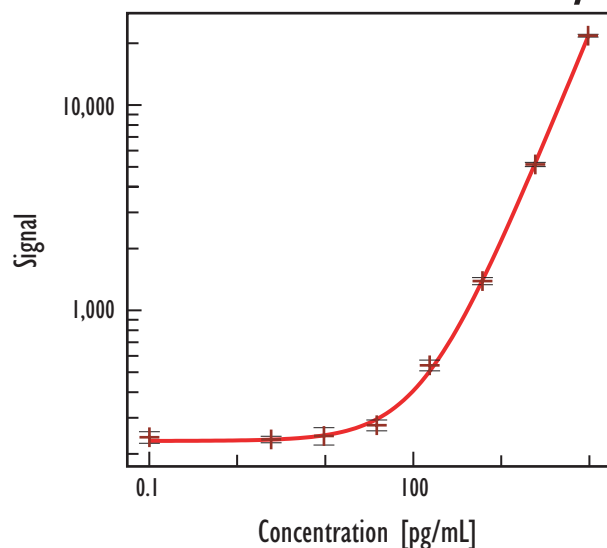
Rat IL-1 $\beta$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	226	13.9
2.4	233	7.0
9.8	253	3.0
39	377	3.1
156	800	3.8
625	2,594	4.2
2,500	10,736	5.9
10,000	47,323	7.3

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	23.5
Serum / Plasma	29.7

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

## Serum / Plasma Assay



Rat IL-1 $\beta$		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	241	6.5
2.4	235	3.7
9.8	245	9.7
39	276	6.1
156	540	6.0
625	1,389	4.0
2,500	5,136	2.4
10,000	21,764	1.6

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	0	28.8 - 123.4
Heparin Plasma (n=10)	1	n/d - 84.7

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.