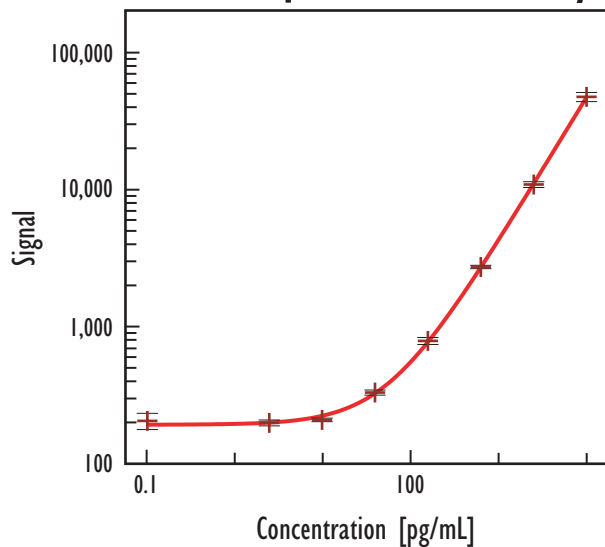


# Rat IL-6 Cytokine Assay (MULTI-SPOT<sup>®</sup> 96-Well 7-Spot Plate)

## Cell Supernatant Assay



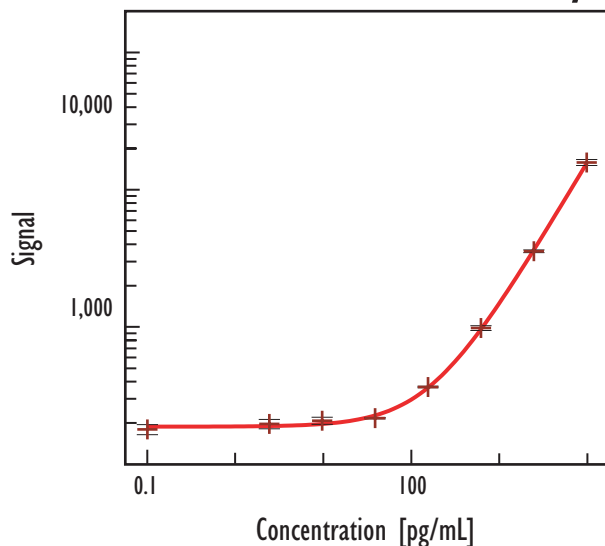
Rat IL-6		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	205	13.6
2.4	198	4.9
9.8	209	3.1
39	330	4.3
156	787	5.9
625	2,727	2.8
2,500	10,902	4.8
10,000	47,559	7.6

## Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	21.3
Serum / Plasma	37.2

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

## Serum / Plasma Assay



Rat IL-6		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	179	8.5
2.4	197	7.8
9.8	208	5.9
39	217	1.1
156	365	1.3
625	982	4.1
2,500	3,560	2.3
10,000	15,791	5.0

## Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	0	49.0 - 194.4
Heparin Plasma (n=10)	5	n/d - 116.8

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.