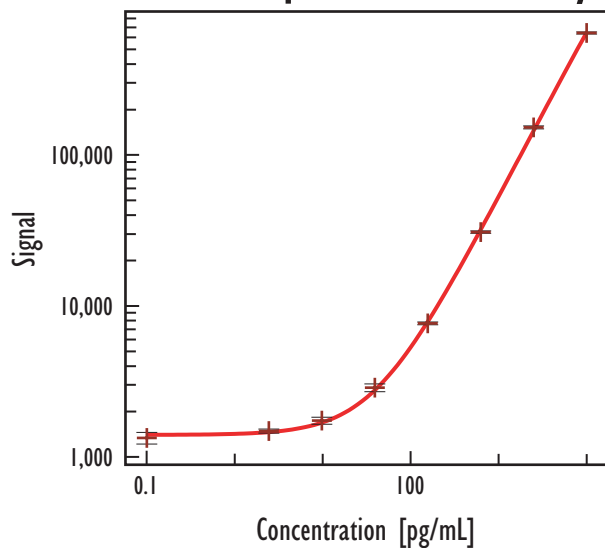


Rat IL-1 α Cytokine Assay (MULTI-SPOT[®] 96-Well 7-Spot Plate)

Cell Supernatant Assay



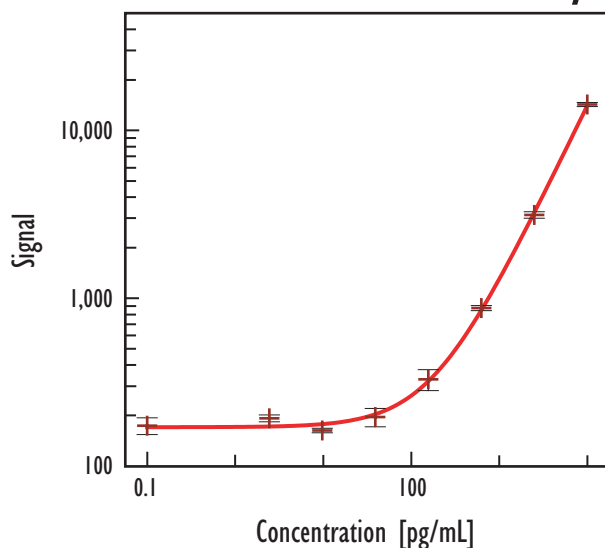
Rat IL-1 α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	1,335	8.8
2.4	1,483	3.1
9.8	1,740	5.3
39	2,875	5.9
156	7,664	2.1
625	30,831	1.8
2,500	152,228	2.3
10,000	642,668	1.6

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	7.5
Serum / Plasma	10.5

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

Serum / Plasma Assay



Rat IL-1 α		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	502	4.6
2.4	554	3.6
9.8	605	7.2
39	905	5.1
156	2,237	1.6
625	7,579	5.1
2,500	33,979	2.0
10,000	161,081	5.7

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=10)	10	n/d
EDTA Plasma (n=10)	5	n/d - 18.8
Heparin Plasma (n=10)	7	n/d - 24.2

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.