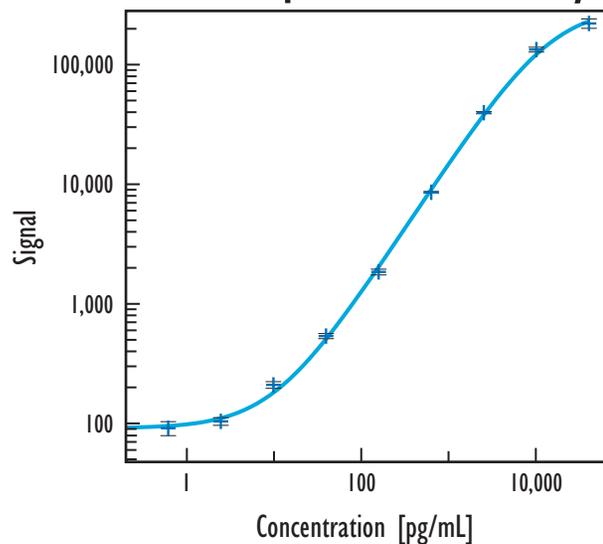


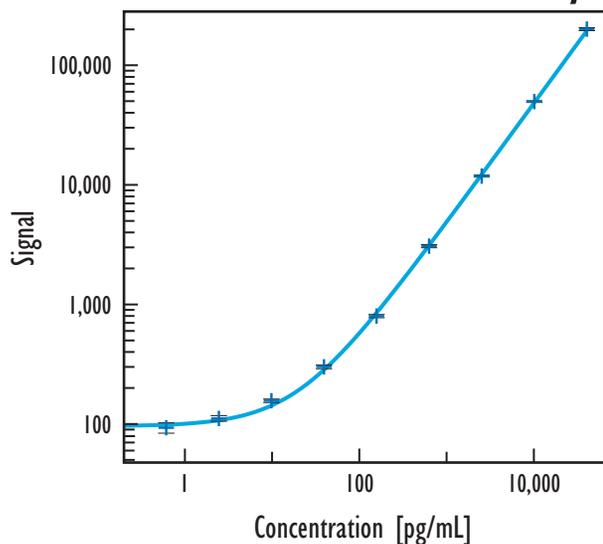
Mouse RANTES Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



Mouse RANTES		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	100	5.9
0.010	92	6.1
0.038	93	5.6
0.15	90	7.8
0.61	92	13.1
2.4	105	7.3
9.8	210	6.2
39	528	4.8
156	1,776	5.4
625	8,033	1.4
2,500	36,210	2.0
10,000	119,509	4.5
40,000	196,067	8.8

Serum / Plasma Assay



Mouse RANTES		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	95	8.6
0.15	93	7.1
0.61	93	10.0
2.4	111	5.4
9.8	156	3.5
39	300	3.7
156	800	3.1
625	3,089	3.1
2,500	11,947	1.9
10,000	50,392	1.7
40,000	204,335	2.9

Detection Limits

Sample Type	Detection Limit (pg/mL)
Cell Supernatant	3.1
Serum / Plasma	5.2

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

Recoveries

Sample Type	Average % Recovery
Cell Supernatant	97
Serum	94
EDTA Plasma	98
Heparin Plasma	100

Spike recoveries were determined in each matrix over a range of spike levels from 39 to 625 pg/mL. Each spike was tested in triplicate.

Endogenous Levels

Sample Type (# of unique samples)	Samples Below Detectable Range	Range
Serum (n=8)	0	77 - 164
EDTA Plasma (n=2)	0	25 - 121
Heparin Plasma (n=8)	0	63 - 140

Endogenous cytokine levels (pg/mL) were determined for different sample types across multiple samples. (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.