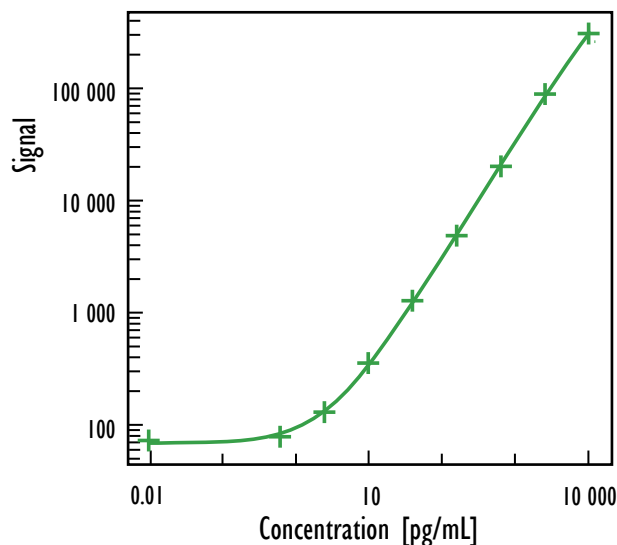


# MULTI-ARRAY<sup>®</sup> Human RANTES Cytokine Assay

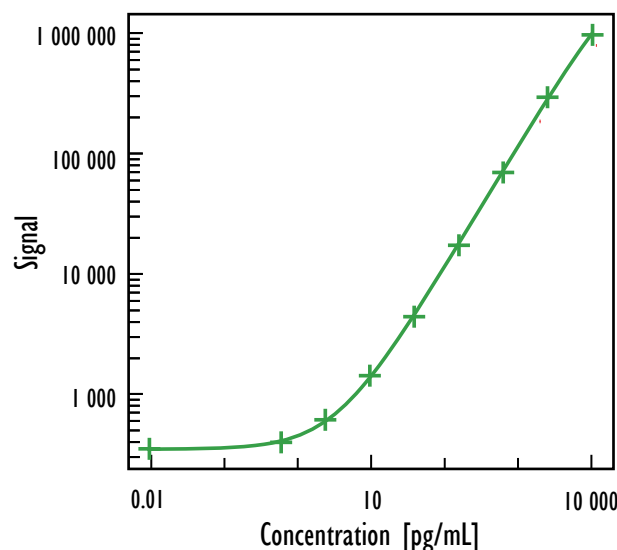
## Tissue Culture Kit



Human RANTES		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	72	7.2
0.61	79	0.7
2.4	132	5.3
9.8	355	5.7
39	1257	5.5
156	4894	3.4
625	20290	4.1
2500	88893	2.4
10000	309801	3.4

Protocol	Detection Limit (pg/mL)
Tissue Culture	0.5

## Ultra-Sensitive Kit



Human RANTES		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	354	2.1
0.61	401	2.3
2.4	616	1.6
9.8	1434	3.5
39	4425	7.7
156	17481	5.3
625	70389	2.4
2500	292137	3.0
10000	969222	0.4

Protocol	Detection Limit (pg/mL)
Serum/Plasma	0.2

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

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K151BFB-1
Tissue Culture	5 plates	K151BFB-2
Tissue Culture	25 plates	K151BFB-4
Ultra-Sensitive	1 plate	K151BFC-1
Ultra-Sensitive	5 plates	K151BFC-2
Ultra-Sensitive	25 plates	K151BFC-4
Base	25 plates	K151BFA-4
Base	250 plates	K151BFA-5

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested. Due to the high endogenous level of Rantes, MSD recommends a dilution of 1:50 fold prior to measuring your sample.

