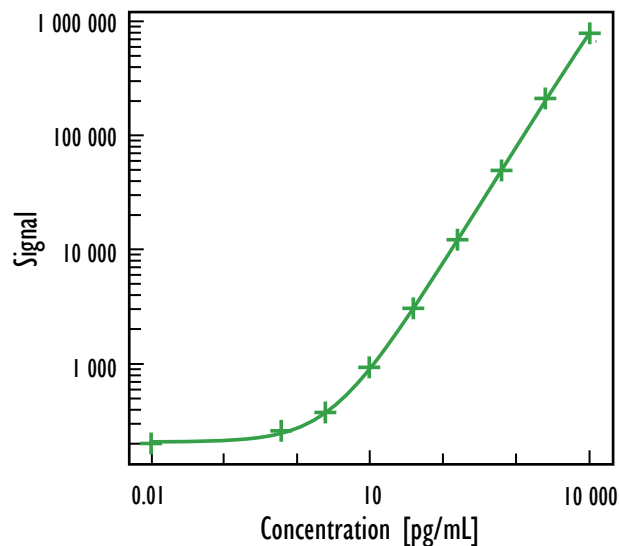


MULTI-ARRAY[®] Human TNF-RII Cytokine Assay

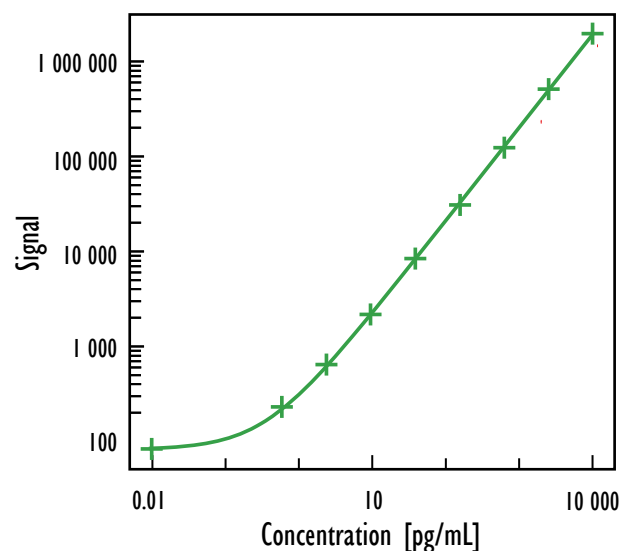
Tissue Culture Kit



Human TNF-RII		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	200	9.0
0.61	255	12.0
2.4	372	2.7
9.8	930	1.2
39	3006	3.3
156	12046	0.7
625	49515	2.2
2500	211608	1.9
10000	790407	1.8

Protocol	Detection Limit (pg/mL)
Tissue Culture	0.5

Ultra-Sensitive Kit



Human TNF-RII		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	83	11.9
0.61	231	3.1
2.4	634	1.2
9.8	2170	0.9
39	8351	1.8
156	30812	3.0
625	124475	1.2
2500	518657	4.1
10000	1976704	3.5

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	K151BJB-1
Tissue Culture	5 plates	K151BJB-2
Tissue Culture	25 plates	K151BJB-4
Ultra-Sensitive	1 plate	K151BJC-1
Ultra-Sensitive	5 plates	K151BJC-2
Ultra-Sensitive	25 plates	K151BJC-4
Base	25 plates	K151BJA-4
Base	250 plates	K151BJA-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 TNFRII, MSD recommends a dilution of 1:10 fold prior to measuring your sample.