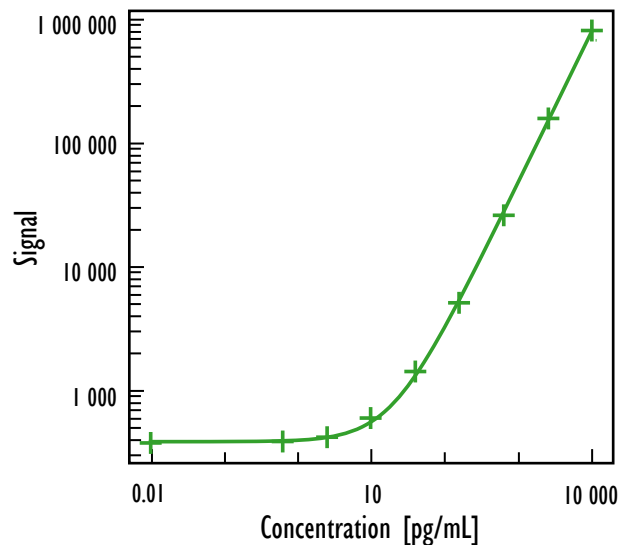


MULTI-ARRAY[®] Human TNF-RI Cytokine Assay

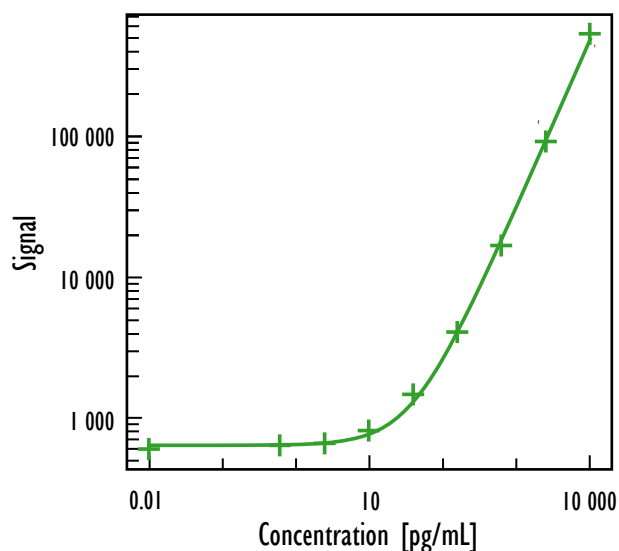
Tissue Culture Kit



Human TNF-RI		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	375	3.2
0.61	390	6.8
2.4	417	6.2
9.8	606	4.1
39	1408	2.1
156	5082	3.8
625	26075	1.3
2500	158100	3.3
10000	825486	3.5

Protocol	Detection Limit (pg/mL)
Tissue Culture	3.2

Ultra-Sensitive Kit



Human TNF-RI		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	606	3.6
0.61	639	5.7
2.4	660	5.8
9.8	814	6.3
39	1503	3.5
156	4095	2.5
625	16566	3.2
2500	92775	3.8
10000	529947	7.1

Protocol	Detection Limit (pg/mL)
Serum/Plasma	8.6

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

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K151BIB-1
Tissue Culture	5 plates	K151BIB-2
Tissue Culture	25 plates	K151BIB-4
Ultra-Sensitive	1 plate	K151BIC-1
Ultra-Sensitive	5 plates	K151BIC-2
Ultra-Sensitive	25 plates	K151BIC-4
Base	25 plates	K151BIA-4
Base	250 plates	K151BIA-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 TNFRI, MSD recommends a dilution of 1:10 fold prior to measuring your sample.