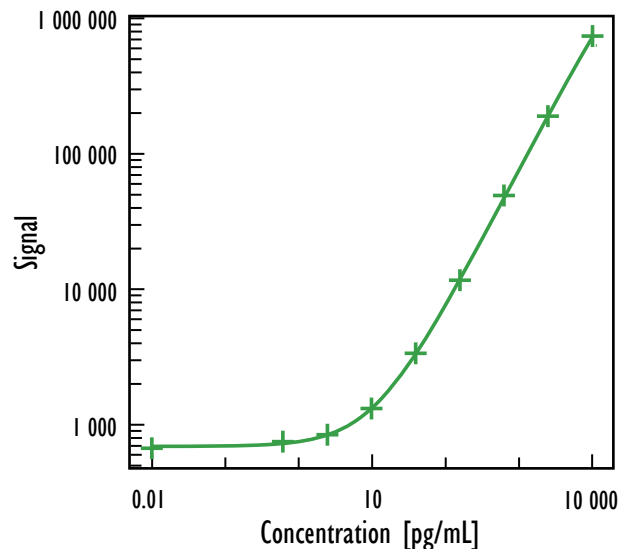


MULTI-ARRAY[®] Human IL-6R Cytokine Assay

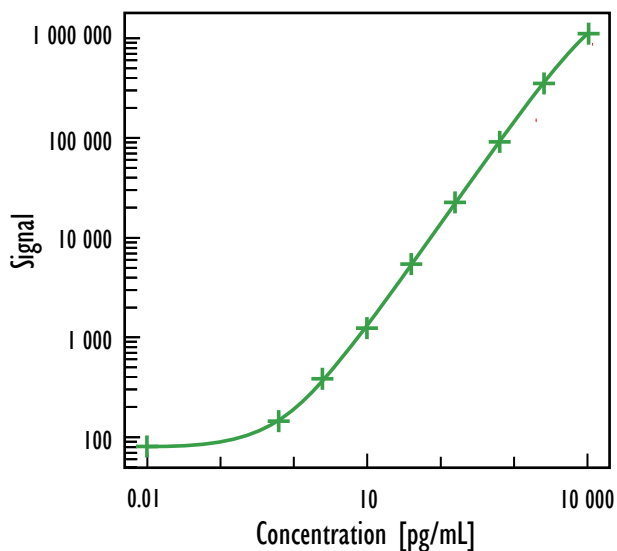
Tissue Culture Kit



Human IL-6R		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	677	5.0
0.61	751	7.2
2.4	847	2.4
9.8	1322	3.5
39	3386	2.5
156	11675	3.9
625	49534	7.0
2500	190802	2.4
10000	737026	4.6

Protocol	Detection Limit (pg/mL)
Tissue Culture	1.4

Ultra-Sensitive Kit



Human IL-6R		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	82	5.2
0.61	142	8.9
2.4	385	1.9
9.8	1252	5.4
39	5399	5.2
156	22620	4.2
625	91925	4.7
2500	352237	4.5
10000	1129331	2.8

Protocol	Detection Limit (pg/mL)
Serum/Plasma	0.1

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

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K151ALB-1
Tissue Culture	5 plates	K151ALB-2
Tissue Culture	25 plates	K151ALB-4
Ultra-Sensitive	1 plate	K151ALC-1
Ultra-Sensitive	5 plates	K151ALC-2
Ultra-Sensitive	25 plates	K151ALC-4
Base	25 plates	K151ALA-4
Base	250 plates	K151ALA-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 IL-6R, MSD recommends a dilution of 1:50 fold prior to measuring your sample.