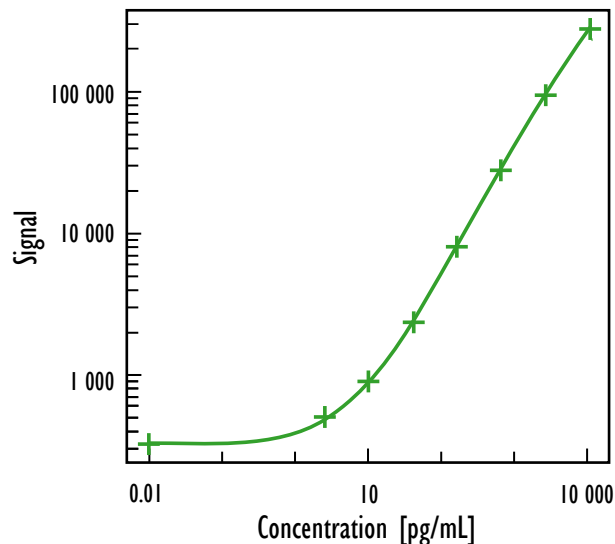


MULTI-ARRAY[®] Human IFN- α 2a Cytokine Assay

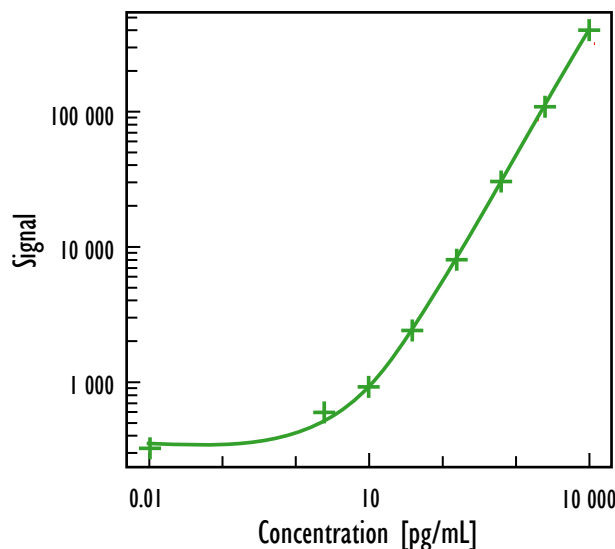
Tissue Culture Kit



Human IFN- α 2a		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	327	1.6
2.4	499	2.2
9.8	887	2.2
39	2355	2.2
156	8038	2.9
625	27856	1.6
2500	94539	1.6
10000	277521	1.7

Protocol	Detection Limit (pg/mL)
Tissue Culture	1.4

Ultra-Sensitive Kit



Human IFN- α 2a		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	324	4.5
2.4	589	4.5
9.8	921	1.7
39	2383	3.4
156	7993	3.4
625	30198	4.5
2500	107315	2.6
10000	396924	2.9

Protocol	Detection Limit (pg/mL)
Serum/Plasma	0.7

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

Kit Model	Kit Size	SI2400	SI6000
Tissue Culture	1 plate	K151ACB-1	K111ACB-1
Tissue Culture	5 plates	K151ACB-2	K111ACB-2
Tissue Culture	25 plates	K151ACB-4	K111ACB-4
Ultra-Sensitive	1 plate	K151ACC-1	K111ACC-1
Ultra-Sensitive	5 plates	K151ACC-2	K111ACC-2
Ultra-Sensitive	25 plates	K151ACC-4	K111ACC-4
Base	25 plates	K151ACA-4	K111ACA-4
Base	250 plates	K151ACA-5	K111ACA-5

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested.