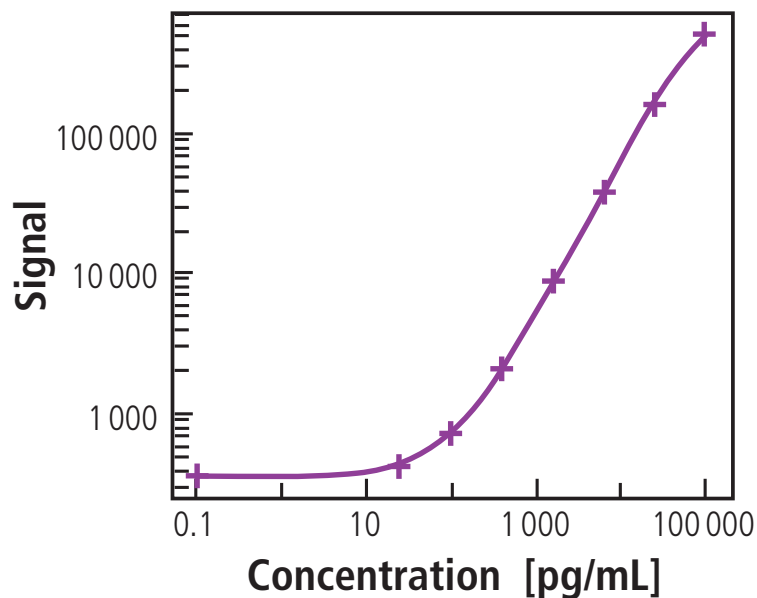


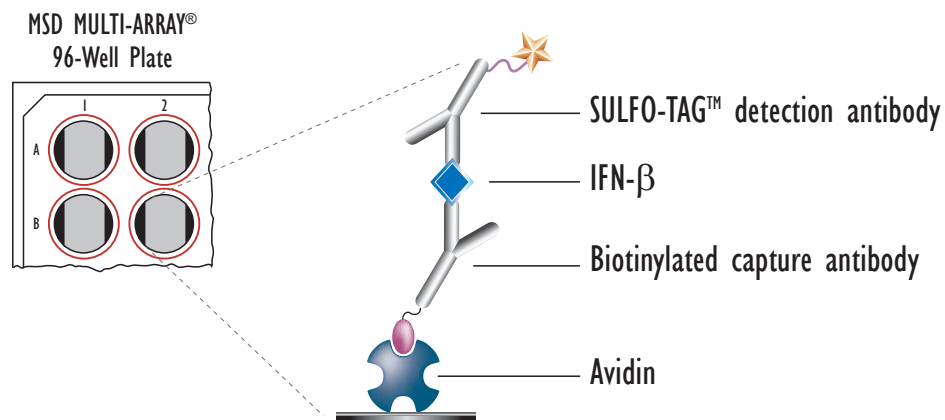
MULTI-ARRAY[®] IFN- β Cytokine Assay

Detection of Human Interferon-beta



Conc (pg/mL)	IFN- β	
	Signal	%CV
0	365	6.5
24	425	2.3
98	711	4.1
391	2107	3.2
1563	8559	1.7
6250	37019	0.3
25000	159748	1.9
100000	486908	1.4

Standard curve data is from a representative experiment



INF- β LLOD	34 (pg/mL)
-------------------	------------

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background

Kit Size	SI2400	SI6000
1 plate	K15IADB-1	K11IADB-1
5 plates	K15IADB-2	K11IADB-2
20 plates	K15IADB-4	K11IADB-4
20 plates (Base)	K15IADA-4	K11IADA-4