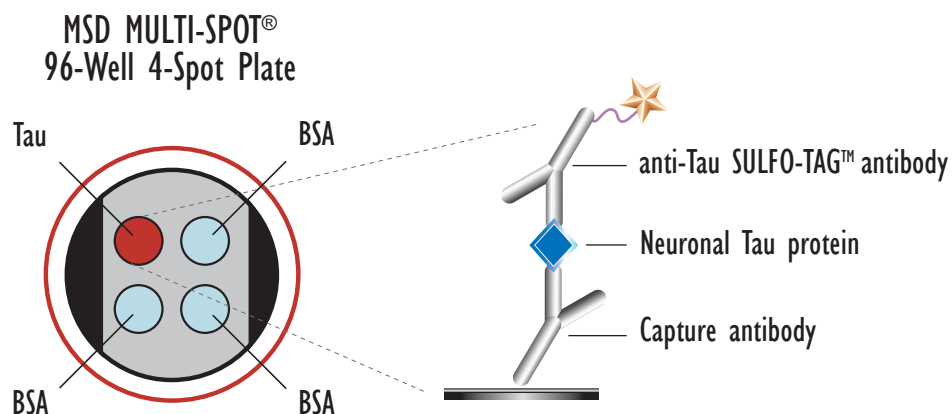
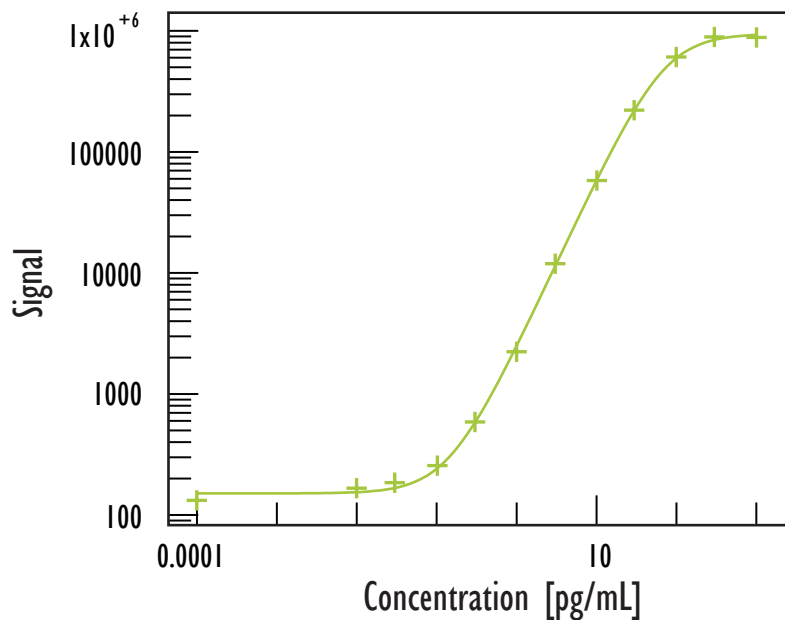


MULTI-ARRAY® Tau Assay

Detection of Tau



ng/mL Tau	Total Tau			S/B
	Average	StdDev	%CV	
0	130	14	11	1
0.01	167	13	8	1
0.03	186	18	10	1
0.1	256	46	18	2
0.3	593	103	17	5
1	2,263	193	9	17
3	11,878	1,112	9	91
10	58,019	5,946	10	445
30	224,690	13,168	6	1,725
100	603,313	30,638	5	4,632
300	878,952	29,606	3	6,748
1,000	889,226	27,203	3	6,827

Total Tau Limit of Detection **35 pg/mL**

Purified neuronal tau was added to MSD MULTI-SPOT 4 Spot 96-well plates coated with anti-total-Tau antibodies on one of the four spatially distinct electrodes per well. Total Tau was detected with anti-total-Tau antibody labeled with MSD SULFO-TAG reagent.

Kit Size	SI2400	SI6000
1 plate	K151DSD-1	K111DSD-1
5 plates	K151DSD-2	K111DSD-2
20 plates	K151DSD-3	K111DSD-3
20 plates (Base)	K151DSA-3	K111DSA-3