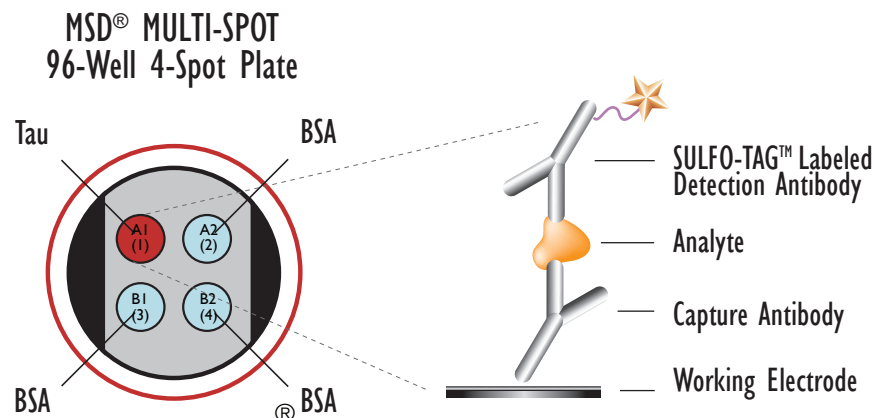
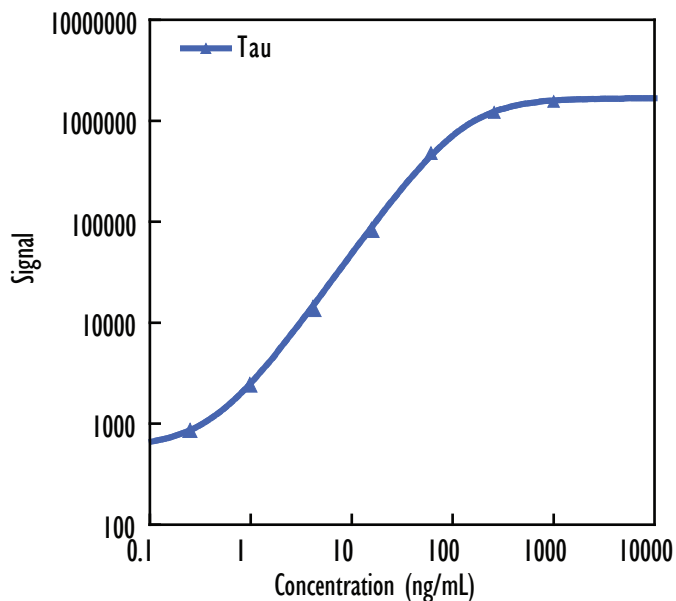


MULTI-ARRAY[®] Mouse Total Tau Assay

Detection of Tau



Tau441 Calibrator (phosphorylated) was added to MSD MULTI-SPOT[®] 4-spot, 96-well plates coated with anti-tau antibody on one of the four spatially distinct electrodes per well. Total tau was detected with anti-tau antibody conjugated with MSD SULFO-TAG reagent.

Tau (ng/mL)	Total Tau	
	Average Signal	%CV
0	483	4.4
0.24	852	1.8
0.98	2440	5.4
3.9	14 020	5.5
16	88 146	2.4
63	465 445	4.4
250	1 256 103	3.7
1000	1 561 670	2.0

Kit Size	Catalog Number
1 plate	K151DSD-1
5 plates	K151DSD-2
20 plates	K151DSD-3
20 plates (Base)	K151DSA-3

Total Tau Limit of Detection	0.076 ng/mL
-------------------------------------	--------------------