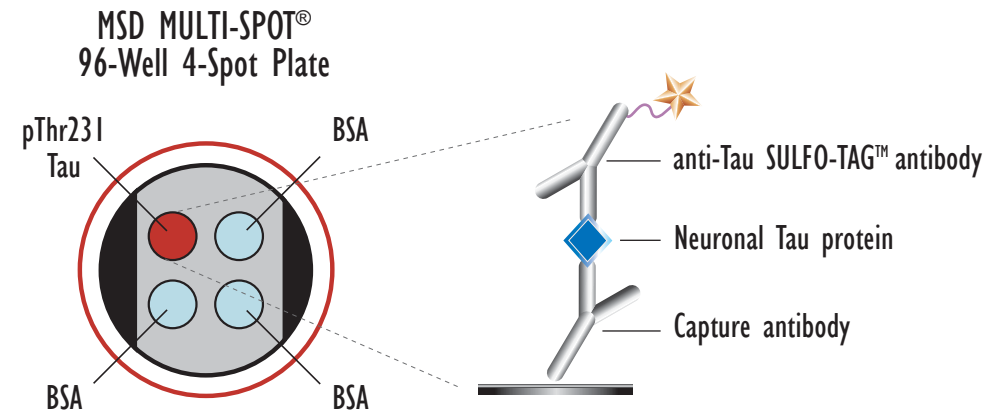
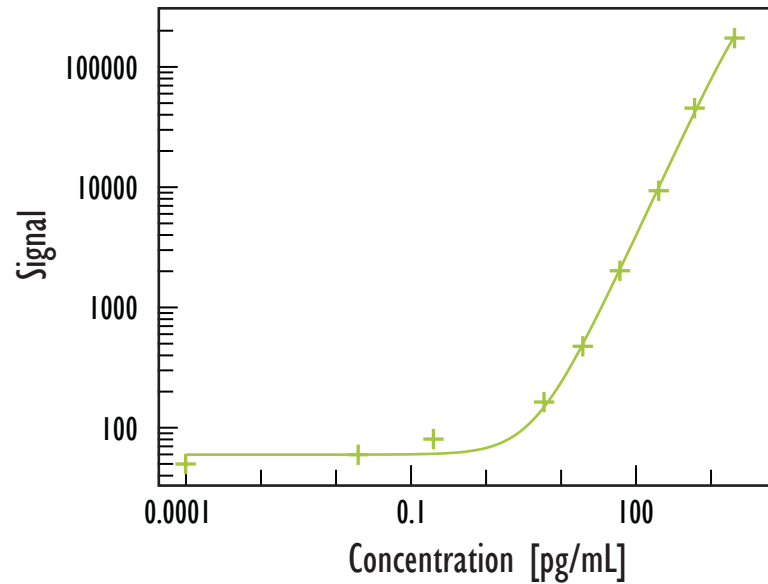


MULTI-ARRAY[®] Phospho- (Thr231) Tau Assay

Detection of Phospho-Tau



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

* Phosphorylated tau is quantified in units because the percentage of the total tau that is phosphorylated may vary with different preparations of the standard.

unit/mL pTau231	Phospho-Tau (Thr231)			S/B
	Average	StdDev	%CV	
0	51	9	17	1
0.02	59	12	20	1
0.2	81	6	7	2
6	167	25	15	3
20	481	27	6	9
60	2,028	93	5	40
200	9,452	517	5	185
600	45,925	2,924	6	900
2000	172,827	6,292	4	3,389

Phospho-Tau (Thr231) Limit of Detection	2.0 units/mL
--------------------------------------------	--------------

Kit Size	SI2400	SI6000
1 plate	K151DRD-1	K111DRD-1
5 plates	K151DRD-2	K111DRD-2
20 plates	K151DRD-3	K111DRD-3
20 plates (Base)	K151DRA-3	K111DRA-3