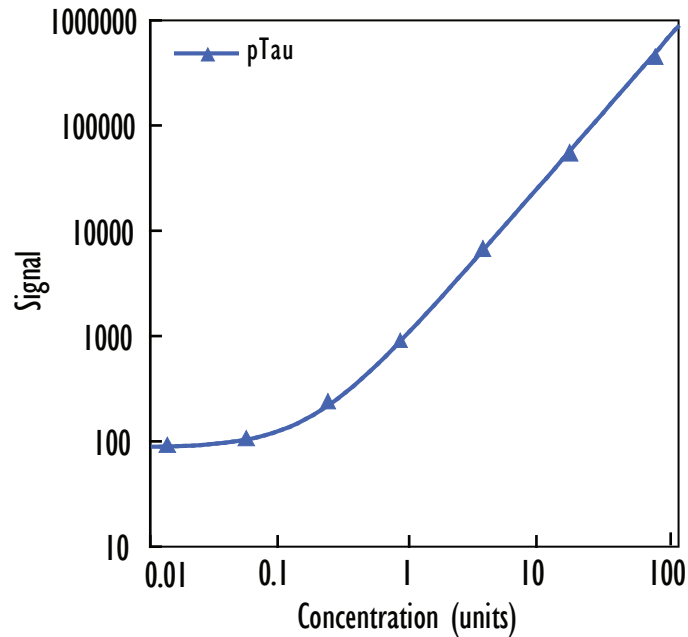


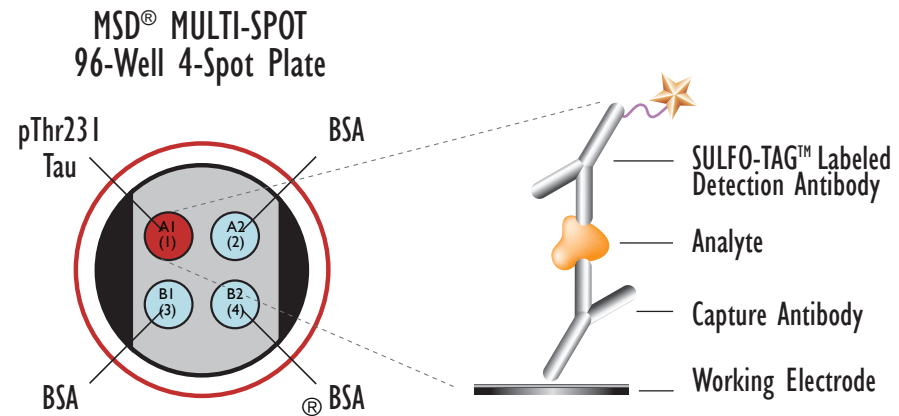
# MULTI-ARRAY<sup>®</sup> Phospho-Tau (Thr231) Assay

## Detection of Phosphorylated Tau



pTau231 (units/well)	Average Signal	%CV
0	78	7.3
0.012	88	9.1
0.049	106	5.2
0.20	217	5.0
0.78	845	4.3
3.1	5384	10.8
13	38 996	4.0
50	326 001	3.5

Phospho-Tau (Thr231) Limit of Detection	0.049 units*
--	--------------



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

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

Kit Size	Catalog Number
1 plate	K151DRD-1
5 plates	K151DRD-2
20 plates	K151DRD-3
20 plates (Base)	K151DRA-3