

AlphaLISA® SureFire® Ultra™

# Human p-Tau (Ser396) Detection Kit

**Product number:** ALSU-PTAU-B500, ALSU-PTAU-B10K,  
ALSU-PTAU-B50K, ALSU-PTAU-B-HV



**Kit specificity:**

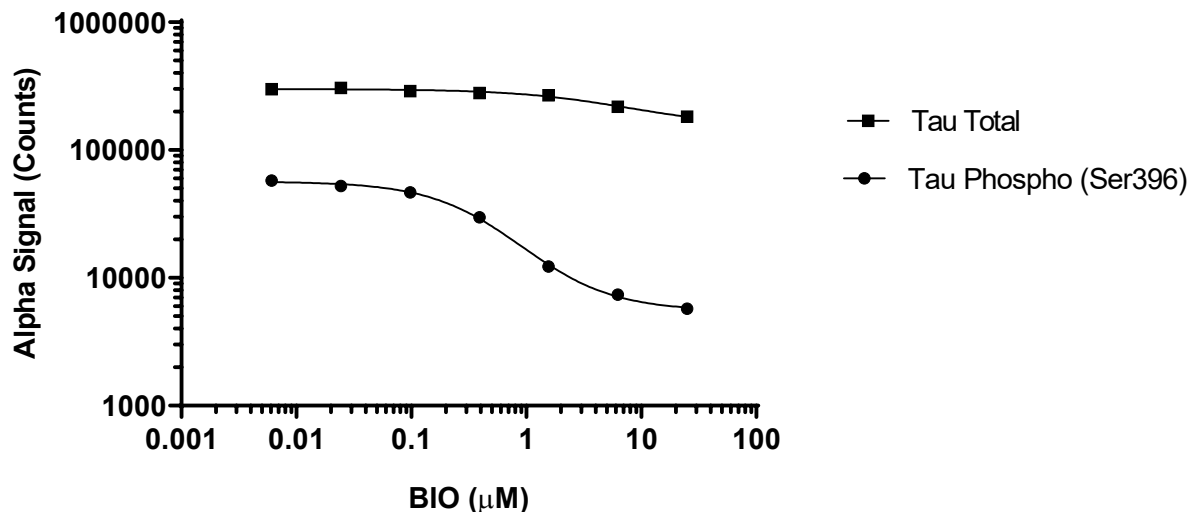
This assay kit contains antibodies which recognize the phospho-Ser396 epitope and a distal epitope on Tau. The protein detected by this kit corresponds to UniProt ID P10636. Tau is also known as Microtubule-associated protein tau. These antibodies recognize Tau of human origin. Other species should be tested on a case-by-case basis.

**Control lysate information:**

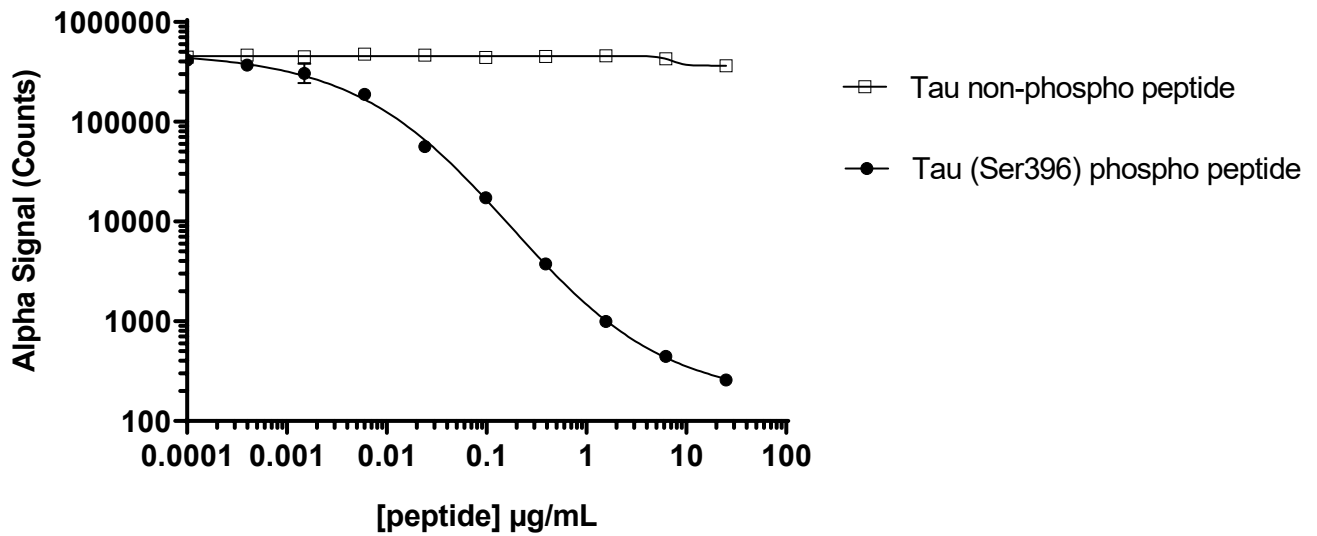
Positive Control Lysate: Prepared from SH-SY5Y cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 100 nM Okadaic Acid for 2.5 hours and lysed with 4 mL of Lysis Buffer.

**Representative data**

Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 20K cells/well in a 96 well plate and incubated for 72 hours. Cells were treated with BIO (GSK-3β inhibitor) prepared in serum free medium, at the indicated concentrations for 3 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser396) and Total Tau using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Phospho (Ser396) or non-phospho Tau peptide were serially diluted into a fixed concentration of SH-SY5Y positive control lysate (25% of stock lysate, equivalent to approx. 12,500 cells/datapoint). Lysates were assayed using the Phospho Tau (Ser396) SureFire Ultra assay kit.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"

