

Instruction Manual

Catalog Number	PK-AB577-3549
Description	Tau, a microtubule-binding protein which serves to stabilize microtubules in growing axons, is found to be hyperphosphorylated in paired helical filaments (PHF), the major fibrous component of neurofibrillary lesions associated with Alzheimer's disease. Hyperphosphorylation of Tau is thought to be the critical event leading to the assembly of PHF. Six Tau protein isoforms have been identified, all of which are phosphorylated by GSK3. This presents the possibility that miscues in GSK3 signaling contribute to the onset of Alzheimer's disease.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 395 of rat Tau.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, chicken, bovine
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-TAU polyclonal antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 0.5% BSA, and 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes mainly the truncated Tau (~30 kDa) in sample from human, mouse and rat origins. Other isoforms of Tau at 35 kDa, 60 kDa and 80 kDa can also be recognized. Reactivity to other species has not been tested. Blocking peptide (Cat. No. PK-AB577-3549P) is available separately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.