

Anti-phospho-Tau (Ser214)

Rabbit Polyclonal Antibody

Catalog # T08-365GR

Lot # O2206-5

Cited Applications

WB

Suggested Dilutions:

WB: 1:500~1:1000

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes endogenous levels of Tau only when phosphorylated at serine 214

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antiserum was produced against synthesized phosphopeptide derived from human Tau around the phosphorylation site of serine 235 (P-K-SP-P-S)

Formulation

PBS + 0.02% sodium azide

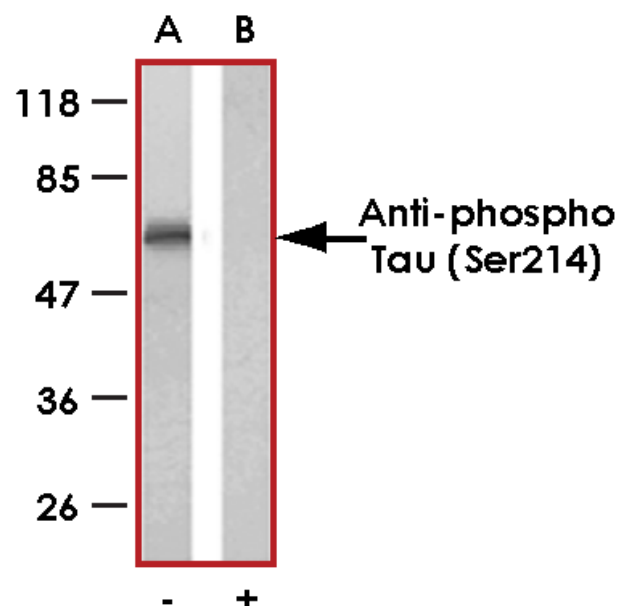
Stability

1yr at -20°C from date of shipment

References

1. Smet C, et al: Regulation of Pin1 peptidyl-prolyl cis/trans isomerase activity by its WW binding module on a multi-phosphorylated peptide of Tau protein. FEBS Lett. 2005 Aug 1;579(19):4159-64.
2. Puig B, et al: Individual and regional variations of phospho-tau species in progressive supranuclear palsy. Acta Neuropathol. 2005 Sep;110(3):261-8.
3. Götz J, et al: Formation of neurofibrillary tangles in P301I tau transgenic mice induced by Abeta 42 fibrils. Science. 2001 Aug 24;293(5534):1491-5.
4. Illenberger S, et al: The endogenous and cell cycle-dependent phosphorylation of tau protein in living cells: implications for Alzheimer's disease. Mol Biol Cell. 1998 Jun;9(6):1495-512.

Sample Data



Western blot analysis of extract from mouse brain tissue using Anti-phospho-Tau (Ser214) antibody

Anti-phospho-Tau (Ser214)

Rabbit Polyclonal Antibody

Catalog Number

T08-365GR

Specific Lot Number

O2206-5

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.