

Anti-phospho-Tau (Ser404)

Rabbit Polyclonal Antibody

Catalog # T08-365HR

Lot # O2206-7

Cited Applications

WB

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 404

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 404 (D-T-SP-P-R)

Formulation

PBS + 0.02% sodium azide

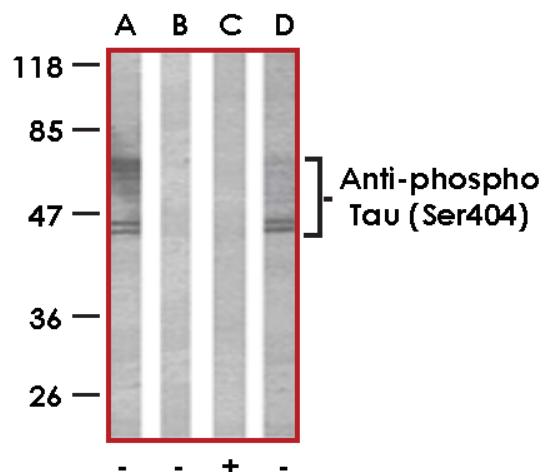
Stability

1yr at -20°C from date of shipment

References

1. Li G, et al: Casein kinase 1 delta phosphorylates tau and disrupts its binding to microtubules. J Biol Chem. 2004 Apr 16;279(16):15938-45.
2. Noble W, et al: Cdk5 is a key factor in tau aggregation and tangle formation in vivo. Neuron. 2003 May 22;38(4):555-65.
3. Giasson BI, et al: The environmental toxin arsenite induces tau hyperphosphorylation. Biochemistry. 2002 Dec 24;41(51):15376-87.
4. Lee G, et al: The microtubule binding domain of tau protein. Neuron. 1989 Jun;2(6):1615-24.
5. Andreadis A et al: Structure and novel exons of the human tau gene. Biochemistry. 1992 Nov 3;31(43):10626-33.

Sample Data



Western blot analysis of extract from mouse brain tissue using a pan-Tau antibody (Lanes A and B) and Anti-phospho-Tau (Ser404) antibody (Lanes C and D)

Anti-phospho-Tau (Ser404)

Rabbit Polyclonal Antibody

Catalog Number

T08-365HR

Specific Lot Number

O2206-7

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.