Phospho-Tau (Ser202, Thr205) recombinant monoclonal antibody (0016)

The anti-Phospho-Tau (Ser202, Thr205) antibody detects phosphorylated Tau protein at serine 202 and threonine 205.

Tau is a microtubule-associated protein that plays a critical role in stabilizing microtubules and supporting axonal transport in neurons. Aberrant phosphorylation of Tau at specific sites, including Ser202 and Thr205, disrupts its normal function and promotes the formation of neurofibrillary tangles (NFTs), a hallmark of Alzheimer's disease and other tauopathies. Phosphorylated Tau at these residues is associated with neuronal dysfunction and disease progression. Hyperphosphorylation of Tau is a key pathological feature in several neurodegenerative diseases, including Alzheimer's disease, frontotemporal dementia, and progressive supranuclear palsy. Detection of Phospho-Tau (Ser202, Thr205) is widely used in research to study disease mechanisms, monitor Tau aggregation, and evaluate potential therapeutic interventions targeting Tau phosphorylation.

This antibody is covered by our Worry-Free Guarantee.

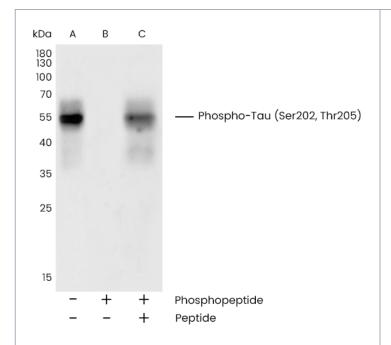
Ordering Information

Order Online »

ENZ-ABS1049-0100 100μl

Manuals, SDS & CofA

View Online »



Western blot analysis of extracts from mouse brain (line A); treated with antigen-specific phosphopeptide (line B) or antigen-specific peptide (line C) using Phospho-Tau (Ser202, Thr205) rabbit monoclonal Antibody (ENZ-ABS1049) at 1:2000 dilution.

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Clone 0016

Formulation Liquid. In 10 mM sodium HEPES, 150 mM NaCl, 100

μg/mL BSA, 50% glycerol, pH 7.5. Preservative-free.

Host Rabbit

Immunogen Synthetic phosphopeptide corresponding to residues

around (Ser202, Thr205) of the Human Phospho-Tau.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/ConditionsWestern Blot (1:2'000-1:10'000)

Suggested dilutions/conditions may not be available for all

applications.

Optimal conditions must be determined individually for

each application.

Species Reactivity Human, Mouse, Rat

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com