

Anti-Phosphothreonine (Anti-pT), Biotin Conjugated

Rabbit Polyclonal Antibody

Catalog # T95-65BR

Lot # E360-6

Cited Applications

For WB (1:1000), ELISA (1:1000)

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

Specificity

Recognizes peptides and proteins phosphorylated on threonine residues.

Cross Reactivity

Pan-specific antibody. No cross-reaction detected with phosphotyrosine peptides and proteins.

Host/Isotype/Clone#

Rabbit

Immunogen

Phosphothreonine peptide-KLH conjugates

Conjugation

Acetylation of primary amine with active ester of biotin (NHS-activated d-biotin).

Formulation

PBS, 50% glycerol, pH7.

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

Scientific Background

Post-translational modification of proteins and peptides is a robust way to regulate function of existing proteins or peptides. Protein phosphorylation represents one of the most abundant and important post-translational modifications of proteins. Phosphorylation on Threonine residues in proteins is performed by a variety of Threonine specific protein kinases in the cell which are involved in a wide variety of signaling networks. The Phospho-Threonine antibody conjugated to Biotin detects phosphorylation on Threonine residues in proteins and peptides.

Anti-Phosphothreonine (Anti-pT), Biotin Conjugated

Rabbit Polyclonal Antibody

Catalog Number

T95-65BR

Specific Lot Number

E360-6

Purification

Affinity chromatography

Concentration

0.25 µg/µL

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.