

Anti-Acetylated Lysine, Biotin Conjugated

Rabbit Polyclonal Antibody

Catalog # L95-66BR

Lot # E372-3

Cited Applications

For WB, ELISA, IP, IHC, IF

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

Specificity

Recognizes acetylated lysine residues in proteins and peptide, including acetylated histone, acetylated BSA and acetylated MBP.

Cross Reactivity

Pan-specific antibody. No cross-reactions detected to the non-acetylated proteins and peptides.

Host/Isotype/Clone#

Rabbit

Immunogen

Acetylated lysine-KLH conjugates

Conjugation

Acylation with NHS-biotin; the biotin : antibody molar ratio was 10:1.

Formulation

PBS, 50% glycerol, pH 7.

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. Acetylation on lysine residues is one example of post-translational modification and is performed by a variety of acetyl transferases in the cell. Histone acetyltransferases (HATs) are universal regulators of eukaryotic transcriptional activity and emerging therapeutic targets for human diseases (1). The Acetylated Lysine Antibody detects acetylation on lysine residues of proteins and peptides.

References

1. Mischerikow, N. et al: Targeted large-scale analysis of protein acetylation. *Proteomics*. 2011 Feb;11(4):571-89.

Sample Data



A B

Representative western blot of the acetylated protein profile with Anti-Acetylated Lysine, Biotin Conjugated (1:1000) using TSA-treated (A) and non-treated (B) human melanoma cells (MMRU) lysate.

Anti-Acetylated Lysine, Biotin Conjugated

Rabbit Polyclonal Antibody

Catalog Number

L95-66BR

Specific Lot Number

E372-3

Purification

By affinity chromatography on an acetyl lysine agarose column

Concentration

0.25 µg/µL

Stability

1 yr 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.