

Acetylated Lysine polyclonal antibody

Acetylation and methylation of lysine are important post-translational modifications that regulate numerous protein-protein and protein-DNA interactions. Lysine acetylation and methylation involves the transfer of acetylCoA, or one or more methyl groups, to the e-amino group of lysine by modifying enzymes and cofactors. Histones and transcription factors are the primary targets of lysine acetylation and methylation, with either modification capable of inducing gene silencing or expression due to differential regulation of cofactors. For example, varying degrees of mono-, di-, and tri-methylation or acetylation of histone H3 at lysine residue 9 are known to demark distinct chromatin regions during various states of gene activation (methylation) or repression (acetylation).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

[View Online »](#)

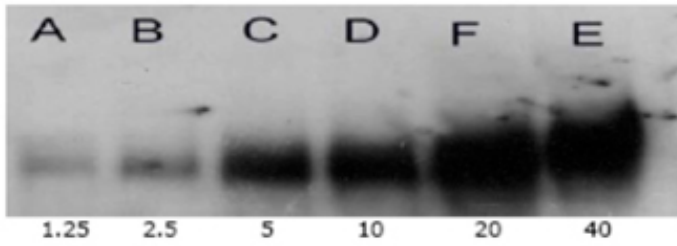
Ordering Information

[Order Online »](#)

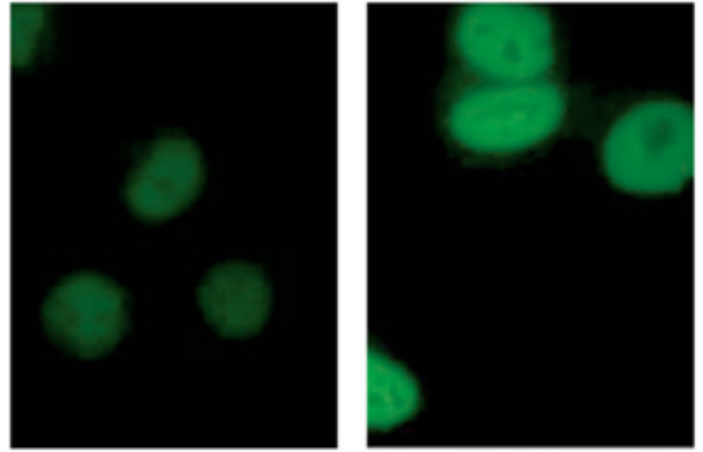
ADI-KAP-TF120-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunoprecipitation analysis of acetylated BSA with Acetylated Lysine (RS1) pAb in ng.



Non-TSA treated

TSA treated

Immunofluorescent analysis of TSA treated (right – 100ng/ml, 24 hours) and non-treated (left) human MMRU melanoma cells using Acetylated Lysine (RS2) pAb.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Ack

Application ELISA, IF, IP, WB

Formulation Liquid. In PBS, 50% glycerol.

Host Rabbit

Immunogen Acetylated KLH.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions ELISA (1:1,000)Immunofluorescence
(1:50)Immunoprecipitation (1:25)Western Blot
(1:250)Suggested dilutions/conditions may not be
available for all applications.Optimal conditions must be
determined individually for each application.

Species Reactivity Species independent

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)