

Methylated Lysine polyclonal antibody (HRP conjugate)

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: 1

[View Online »](#)

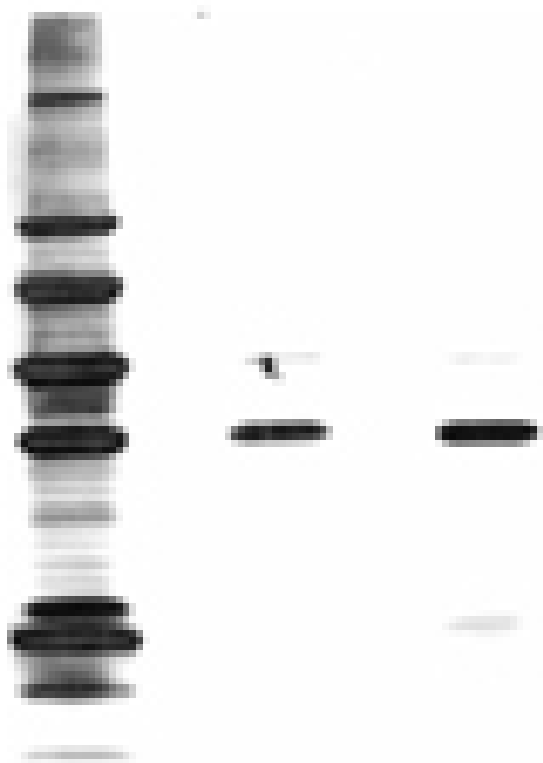
Ordering Information

[Order Online »](#)

ADI-KAP-TF1213-E	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of the methylated protein pattern in whole HeLa cell with Methylated Lysine pAb, HRP conjugate. Left: Signal of methylated proteins; Middle: Signal of the methylated proteins was competitively suppressed by 50µg/ml of methylated BSA; Right: Signal of the methylated proteins was competitively suppressed by 10µg/ml of methylated BSA.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, IHC, WB

Formulation Liquid. In HRP stabilizing buffer with 50% glycerol.

Host Rabbit

Immunogen Methylated KLH.

Purity Detail Immunoaffinity purified.

Recommendation Dilutions/Conditions ELISA (0.1µg/ml)Western Blot (0.2-0.5µg/ml)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-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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