

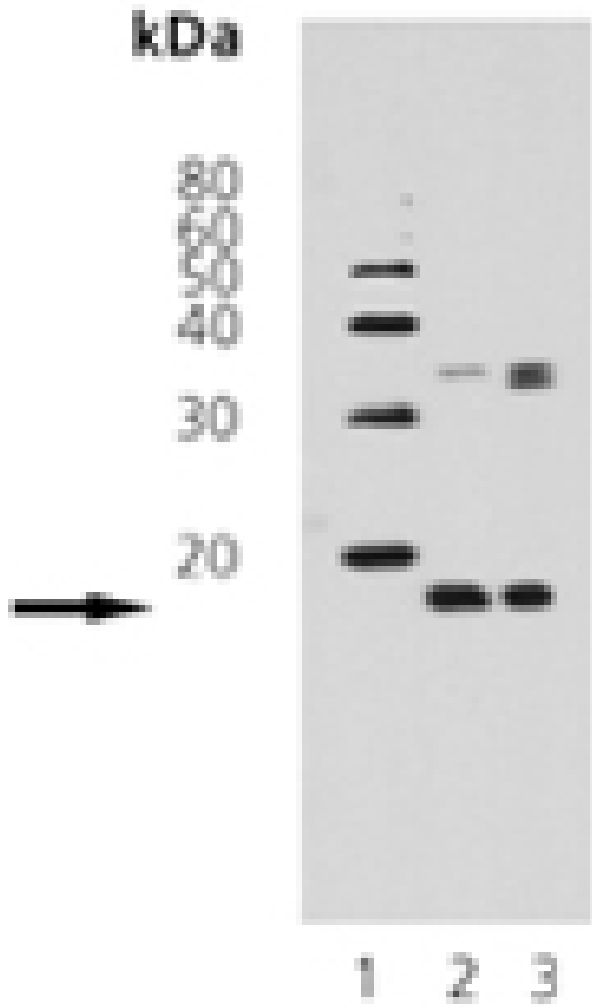
Histone H3 (dimethyl-Lys9) polyclonal antibody

The nucleosome is made up of four core histone proteins (H2A, H2B, H3 and H4) and is the primary building block of chromatin. The core histones form an octamer, which is made up of one H3-H4 tetramer and two H2A-H2B dimers. The N-terminal tail of core histones undergoes different post-transcriptional modifications including acetylation, phosphorylation, methylation and ubiquitination. These modifications occur in response to cell signal stimuli and have a direct effect on gene expression.

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

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of MW marker (1), 3T3 (2), and K-562 (3) cell lysates probed with Histone H3 (dimethyl-Lys9) pAb.

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

Crossreactivity Weakly cross-reacts with rat, bovine and dog Histone H3 (dimethyl-Lys9).

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Host Rabbit

Immunogen Synthetic dimethyl-peptide corresponding to human Histone H3.

Purity Detail Peptide affinity purified.

Source Purified from rabbit serum.

Species Reactivity Human, Mouse

UniProt ID P68431 (Histone H3.1), Q71DI3 (Histone H3.2), P84243 (Histone H3.2)

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)