## 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.

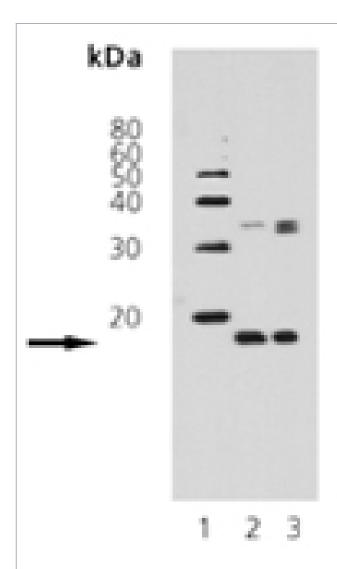
**Ordering Information** 

Order Online »

**ADI-905-778-100** 100μg

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com