

## JMJD2A antibody (mAb)

**Catalog Nos:** 39815, 39816**RRID:** AB\_2793352**Clone:** N154/32**Isotype:** IgG1**Application(s):** WB**Reactivity:** Human, Mouse**Quantities:** 100 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Concentration:** 1 µg/µl**Molecular Weight:** 130 kDa

**Background:** JMJD2A (KDM4A, JHDM3A) is a member of the Jmj-containing (Jumonji) class of histone demethylase proteins that are involved in the regulation of genome function through the removal of methyl groups from histones. JMJD2A specifically demethylates histone H3 dimethyl lysine 9 (H3K9me2), trimethyl lysine 9, (H3K9me3), dimethyl lysine 36 (H3K36me2) and trimethyl lysine 36 (H3K36me3). It also participates in transcriptional repression of ASCL2 and E2F-responsive promoters via the recruitment of histone deacetylases and NCOR1. JMJD2A contains JmjC and JmjN domains, two PHD fingers and two tudor domains.

**Immunogen:** This JMJD2A antibody was raised against a recombinant protein corresponding to amino acids 586-674 of human JMJD2A.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

Applications Validated by Active Motif:

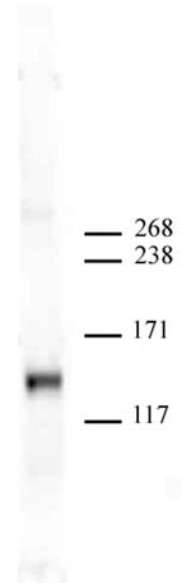
WB\*: 1 - 2 µg/ml dilution

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot. The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is also recommended to aid in detection by Western blot.

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91143.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**JMJD2A antibody (mAb) tested by Western blot.**

HeLa cell nuclear extract (40 µg) probed with JMJD2A antibody at a dilution of 2 µg/ml.