

JMJD2D antibody (pAb)

Catalog Nos: 39247, 39248**RRID:** AB_2793203**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 58 kDa

Background: JMJD2D (KDM4D) is a histone demethylase located in the nucleus that specifically demethylates Lysine 9 of Histone H3, thereby playing a central role in the histone code. JMJD2D (KDM4D) demethylates both di- and trimethylated histone H3 Lys 9, while it has no activity on monomethylated residues. The demethylation reaction of the Lys residue generates formaldehyde and succinate as a by product. In addition, JMJD2D forms a complex with the ligand-bound form of the Androgen Receptor (AR) and, through this interaction, activates AR expression. Overexpression of AR is associated with prostate cancer, suggesting that, via its ability to upregulate AR, JMJD2D (KDM4D) may be involved in carcinogenesis.

Immunogen: This JMJD2D antibody was raised against a peptide corresponding to amino acids 337-359 of human JMJD2D.

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

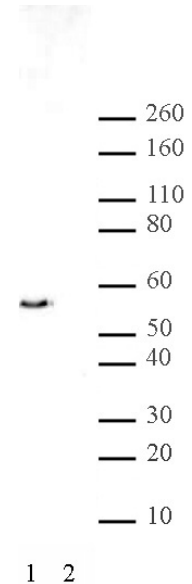
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 dilution

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.

**JMJD2D pAb tested by Western blot.**

Nuclear extract of HeLa cells (20 µg) probed with JMJD2D pAb (1:1,000). Specificity was verified by peptide competition.

Lane 1: No peptide.

Lane 2: 1 µM immunizing peptide.