

JMJD6 antibody (pAb)

Catalog Nos: 61493, 61494**RRID:** AB_2793657**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 52 kDa

Background: JMJD6 (Jumonji Domain-Containing Protein 6) is a dioxygenase that can both act as a histone arginine demethylase and a lysyl-hydroxylase. JMJD6 acts as an arginine demethylase which demethylates histone H3 at 'Arg -2' (H3R2me) and histone H4 at 'Arg-3' (H4R3me), thereby playing a role in histone code. Required for differentiation of multiple organs during embryogenesis and acts as a key regulator of hematopoietic differentiation. May act as a RNA hydroxylase, as suggested by its ability to bind single strand RNA.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human JMJD6.

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:

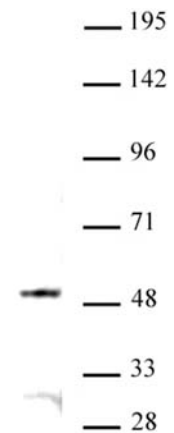
WB*: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*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.

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.

**JMJD6 antibody (pAb) tested by Western blot.**

Nuclear extract of Raji cells (20 µg) probed with JMJD6 antibody (pAb) at a dilution of 1:500.