

KDM4C antibody (pAb)

Catalog Nos: 61687, 61688

RRID: AB_2793735

Isotype: IgG

Application(s): WB

Reactivity: Monkey

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 142 kDa

Background: **KDM4C** (Lysine (K)-Specific Demethylase 4C) is a histone demethylase that specifically demethylates 'Lys-9' and 'Lys-36' residues of histone H3, thereby playing a central role in histone code. Does not demethylate histone H3 'Lys-4', H3 'Lys-27' nor H4 'Lys-20'. Demethylates trimethylated H3 'Lys-9' and H3 'Lys-36' residue, while it has no activity on mono- and dimethylated residues. Demethylation of Lys residue generates formaldehyde and succinate.

Immunogen: This antibody was raised against a peptide within human KDM4C.

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

Application Notes:

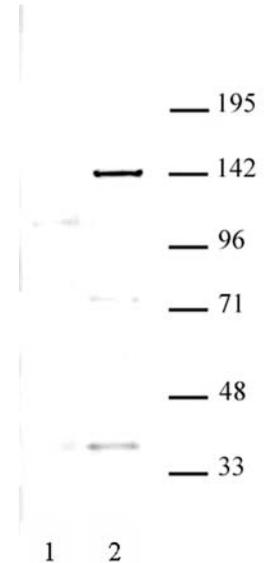
Applications Validated by Active Motif:

WB: 1:500 dilution

For JMJD2C / KDM4C, we also offer AbFlex[®] JMJD2C / KDM4C Recombinant Antibody (rAb). For details, see Catalog No. 91307.

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.



KDM4C antibody (pAb) tested by Western blot.

COS-7 nuclear extract (25 µg per lane) was probed with KDM4C pAb at a 1:500 dilution. Lane 1: No treatment. Lane 2: Cells treated with CoCl₂.