

## KDM4B Polyclonal Antibody

(Catalog # A-2727)

### Background

Histone demethylase that specifically demethylates 'Lys-9' of histone H3, thereby playing a role in histone code. Does not demethylate histone H3 'Lys-4', H3 'Lys-27', H3 'Lys-36' nor H4 'Lys-20'. Only able to demethylate trimethylated H3 'Lys-9', with a weaker activity than KDM4A, KDM4C and KDM4D. Demethylation of Lys residue generates formaldehyde and succinate.

### Description

KDM4B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

O94953

### Purification

Affinity Purified

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 917-1031 of human KDM4B/JMJD2B (NP\_055830.1)

### Storage

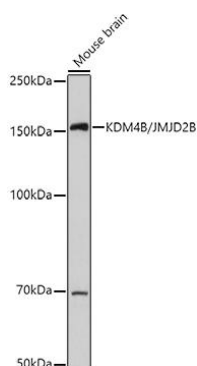
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

JMJD2B, TDRD14B

### Application

WB; Recommended dilution: WB 1:500 - 1:1000



Western blot analysis of extracts of Mouse brain, using KDM4B Polyclonal Antibody at 1:1000 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.