

KDM2B antibody (mAb)

Catalog Nos: 65699, 65999, 65700

RRID: AB_3216371

Clone: K2K842-13

Application(s): WB

Reactivity: Human, Mouse

Quantities: 100 µg, 50 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Isotype: IgG

Molecular Weight: 100 kDa

Background: KDM2B (lysine (K)-specific demethylase 2B), also known as FBXL10 (F-box and leucine-rich repeat protein 10) is a histone demethylase that preferentially demethylates trimethylated lysine 4 (K4me3) and dimethylated lysine 36 (K36me2) of histone H3. KDM2B displays weak or no activity for mono- and trimethylated H3K36. KDM2B preferentially binds the transcribed region of ribosomal RNA and represses the transcription of ribosomal RNA genes which results in inhibition of cell growth and proliferation. KDM2B may also serve as a substrate-recognition component of the SCF (SKP1-CUL1-F-box protein)-type E3 ubiquitin ligase complex.

Immunogen: This KDM2B antibody was raised against a peptide containing human lysine-specific demethylase 2B of isoform a.

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

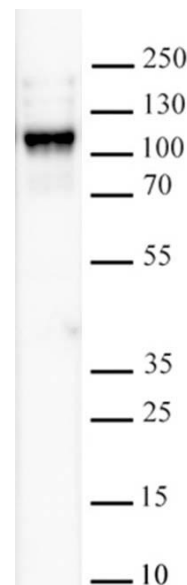
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml 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.



KDM2B antibody (mAb) tested by Western blot.

Detection of KDM2B by Western blot analysis.

20 µg of HEK293 nuclear extract was run on SDS-PAGE and probed with 0.5 µg of KDM2B antibody.