

JARID1B / KDM5B antibody (pAb)

Catalog Nos: 65730, 65930, 65731**RRID:** AB_3216312**Application(s):** ChIP**Reactivity:** Human**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 180 kDa

Background: Lysine (K)-specific demethylase 5B (KDM5B), also known as Jumonji, AT rich interactive domain 1B (JARID1B), is a histone demethylase that specifically demethylates lysine 4 of histone H3 (H3K4), thereby playing a central role in defining the histone code. KDM5B demethylates trimethylated, dimethylated and monomethylated H3K4. It does not demethylate histone H3K9 or H3K27. KDM5B functions as a transcriptional corepressor for FOXG1B and PAX9. KDM5B induces the proliferation of breast cancer cells by repressing tumor suppressor genes such as BRCA1 and HOXA5. In contrast, it may act as a tumor suppressor for melanoma.

Immunogen: This antibody was raised against a human peptide belonging to demethylase JARID1B.

Buffer: Purified IgG in PBS, 30% glycerol, 0.035% Azide. Sodium azide is highly toxic.

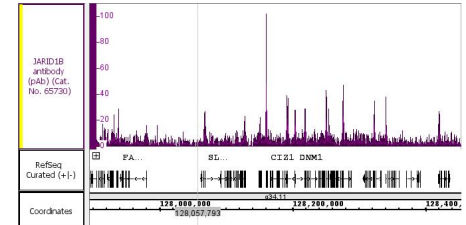
Application Notes:

Applications Validated by Active Motif:

ChIP-Seq: 10 µl each

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.

**JARID1B / KDM5B antibody (pAb) tested by ChIP-Seq.**

ChIP was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 40 µg of chromatin from K562 cells and 10 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify JARID1B / KDM5B binding sites. The image shows binding across a region of chromosome 9. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.