

JARID1C / KDM5C antibody (pAb)

Catalog Nos: 39263, 39264**RRID:** AB_2615032**Isotype:** IgG**Application(s):** WB**Reactivity:** Budding Yeast**Volumes:** 200 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 85 kDa

Background: JARID1C / KDM5C is a JmjC domain family histone demethylase specific for H3K4 (lysine at position 4 of the histone H3 protein). It removes methyl groups specifically added by Set1p methyltransferase.

Immunogen: This antibody was raised against a peptide corresponding to the C-terminus of JARID1C / KDM5C from *S. cerevisiae*.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol, 0.7 mg/ml BSA and 0.035% sodium azide. Sodium azide is highly toxic.

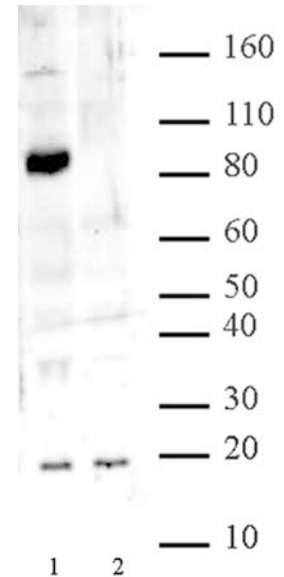
Application Notes:

Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 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.

**JARID1C / KDM5C pAb tested by Western blot**

S. cerevisiae extracts probed with JARID1C / KDM5C pAb (1:1,000 dilution).

Lane 1: wild type strain.

Lane 2: JARID1C / KDM5C knockout strain.