

ALKBH5 antibody (rAb), 100 µl

Catalog No: 91573

Clone: 2K9

Application(s): IF, WB

Reactivity: Human, Mouse, Rat

Volume: 100 µl

Purification: Protein A Chromatography

Host: Rabbit

Isotype: IgG

Molecular Weight: 57 kDa

Background: **ALKBH5** (ABH5 and OFOX1) is a Dioxygenase that demethylates RNA by oxidative demethylation. It is a mammalian demethylase that oxidatively reverses m6A in mRNA in vitro and in vivo. **ALKBH5** is required for spermatogenesis. Two major isoforms for this target are located at a of MW 40-44 kDa and 52 kDa.

Immunogen: This antibody was raised against the ALKBH5 fusion protein.

Buffer: Purified IgG in PBS, pH 7.3, with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

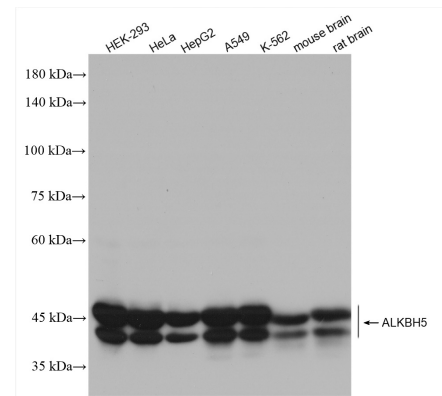
Applications Validated by Active Motif:

WB: 1:5000-1:50000 dilution

IF: 1:250-1:1000 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.



ALKBH5 antibody (rAb) tested by Western blot.

Detection of ALKBH5 by Western blot analysis using various lysates and probed with ALKBH5 antibody at 1:10000 dilution.