

AbFlex® ALKBH5 antibody (rAb)

Catalog Nos: 91441, 91442

Application(s): IF, IP, WB

Reactivity: Human, Mouse, Rat

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Isotype: IgG

Molecular Weight: 44 kDa

Background: AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports.

AbFlex® ALKBH5 antibody was expressed as full-length IgG with rabbit immunoglobulin heavy and light chains in mammalian HEK293 cells. **ALKBH5** (AlkB Homolog 5, RNA Demethylase, also known as OFOXD1, ABH5) specifically demethylates m6A RNA. **ALKBH5** significantly affects mRNA export and RNA metabolism as well as the assembly of mRNA processing factors in nuclear speckles via demethylation activity.

Immunogen: This ALKBH5 antibody was raised against a peptide of human ALKBH5.

Buffer: Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

*IP: 2 µg per IP

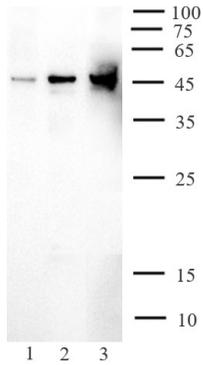
WB: 0.5 - 2 µg/ml dilution

IF: 2 µg per ml

*For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

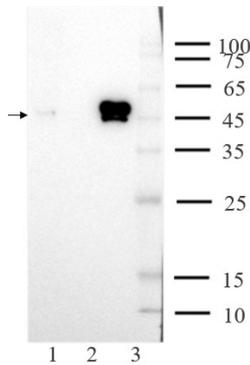
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



AbFlex® ALKBH5 rAb antibody tested by Western blot.

Detection of ALKBH5 by Western blot analysis using 50 µg of HepG2 (Lane 1), A375 (Lane 2) and HEK293 (Lane 3) whole cell extracts and probed with antibody at 0.5 µg/ml.



AbFlex® ALKBH5 rAb antibody tested by immunoprecipitation.

2 µg of antibody was used to immunoprecipitated ALKBH5 from 1 mg of HEK293F whole cell extract, 40% IP product (Lane 3) and 40 µg of Input proteins (Lane 1) were then run on SDS-PAGE and probed with the same antibody at 0.5 µg /ml. Tween-20 at 0.1% was added to blocking buffers to perform Western blot and reduce background. 2 µg of rabbit IgG was used as negative control for IP (Lane 2).

AbFlex® ALKBH5 rAb antibody tested by immunofluorescence.

Staining of HeLa cells with ALKBH5 antibody at 2 µg/mL. Followed by Multi-rAb CoraLite® Plus 594-Goat Anti-Rabbit Recombinant Secondary Antibody (H+L). DAPI at 594 nm

