

FTO antibody (pAb)

Catalog Nos: 61431, 61432**RRID:** AB_2793633**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 60 kDa

Background: FTO (fat mass and obesity associated) is a dioxygenase that repairs alkylated DNA and RNA by oxidative demethylation. This protein has high activity towards single-stranded RNA containing 3-methyluracil and single-stranded DNA containing 3-methylthymine. FTO has low demethylase activity towards single-stranded DNA containing 1-methyladenine or 3-methylcytosine and has no activity towards 1-methylguanine or double-stranded DNA. FTO contributes to the regulation of the global metabolic rate, energy expenditure, energy homeostasis, regulation of body size and body fat accumulation.

FTO has recently been shown to have the highest activity towards demethylation of 6-methyladenine in single-stranded RNA suggesting that FTO is involved in processing of nuclear RNAs and may play an important role in RNA epigenetics.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human FTO.

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

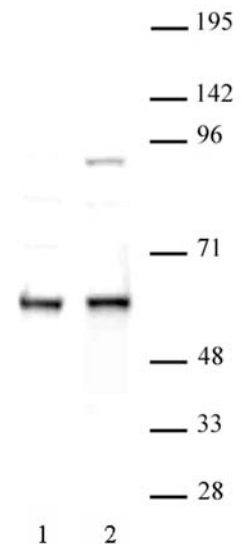
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 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.



FTO antibody (pAb) tested by Western blot. Detection of FTO by Western blot analysis (30 µg per lane). Lane 1: Nuclear extract of Caco-2 cells. Lane 2: Nuclear extract of HEK-293 cells. Both probed with FTO antibody (pAb) at a 1:500 dilution.