

APOBEC2 antibody (pAb)

Catalog Nos: 39985, 39986

RRID: AB_2793418

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 35 kDa

Background: APOBEC2 (Apolipoprotein B mRNA-editing enzyme catalytic subunit 2) is part of a family of enzymes that edit nucleic acids, possibly changing cytosine to uracil. APOBEC2 expression is regulated by TNF- α via NF κ B activation and this suggests a possible role for APOBEC2 in the pathophysiology of hepatic inflammation. APOBEC2 has a low intrinsic cytidine deaminase activity and is also involved in DNA demethylation. APOBEC2 deaminates 5-methylcytosine, followed by base excision repair by Mbd4.

Immunogen: This APOBEC2 antibody was raised against a peptide derived from the N-terminal region of human APOBEC2.

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

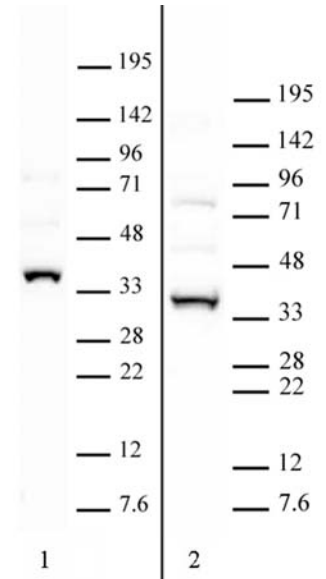
Application Notes:

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.



APOBEC2 antibody (pAb) tested by Western blot.

Nuclear extract (20 μ g) of L6E9 cells (Lane 1) or nuclear extract (20 μ g) of C2C12 cells (Lane 2) probed with APOBEC2 antibody at a 1:2,000 dilution.