

Cxxc4 antibody (pAb)

Catalog Nos: 61679, 61680

RRID: AB_2793732

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 42 kDa

Background: **Cxxc4** (CXXC Finger Protein 4) acts as a negative regulator of the Wnt signaling pathway via its interaction with DVL1 and has been implicated in malignant renal cell carcinoma and colonic villous adenoma. Recently, Cxxc4 has been shown to be involved in the regulation of DNA demethylation. It binds to unmethylated CpGs in DNA and is involved in TET2 downregulation.

Immunogen: This antibody was raised against a peptide within the N-terminal region of mouse Cxxc4.

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

Application Notes:

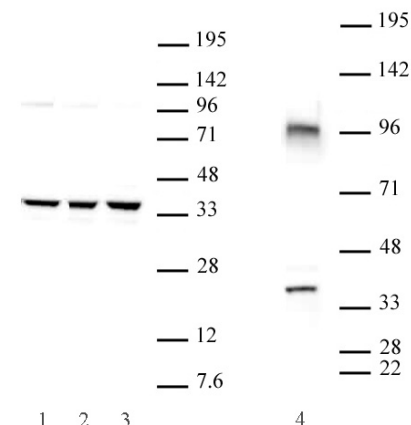
Applications Validated by Active Motif:

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

The addition of 0.1% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. 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.



Cxxc4 antibody (pAb) tested by Western blot.

Detection of Cxxc4 antibody by Western blot analysis.

Lane 1: Mouse ES cytoplasmic extract.

Lane 2: F9 cytoplasmic extract.

Lane 3: P19 cytoplasmic extract.

Lane 4: HEK293 whole cell extract.

30 µg of lysate per lane probed with Cxxc4 antibody at a dilution of 1:1,000.