

C17orf96 antibody (pAb)

Catalog Nos: 61753, 61754

RRID: AB_2793758

Isotype: IgG

Application(s): ChIP-Seq, WB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 kDa

Background: C17orf96 has been shown to be highly expressed in mammalian ES cells and has been implicated as an interacting protein of the PRC2 complex (polycomb repressive complex 2). It may function as a CpG island binding protein and thus a potential regulator of gene transcription.

Immunogen: This antibody was raised against a recombinant protein corresponding to amino acids 59-229 of mouse C17orf96.

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.

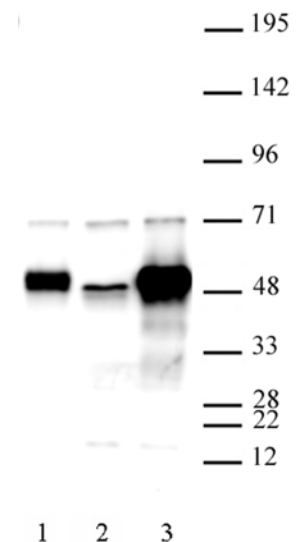
Application Notes:

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.



C17orf96 antibody (pAb) tested by Western blot.

Detection of C17orf96 by Western blot analysis.

Lane 1: Mouse ES cell nuclear extract.

Lane 2: Mouse ES cell nuclear extract from cells with a C17orf96 knockdown.

Lane 3: Mouse ES cell nuclear extract from cells with C17orf96 overexpressed.

All probed with C17orf96 antibody at a dilution of 1:1,000.