

CHD2 antibody (pAb)

Catalog Nos: 39363, 39364

RRID: AB_2793228

Isotype: IgG

Application(s): WB

Reactivity: Human, Wide Range Predicted

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 260 kDa

Background: CHD (Chromo-ATPase/Helicase-DNA-binding) proteins, which are ATP-dependent chromatin remodeling factors implicated in RNA polymerase II transcriptional regulation, are characterized by the presence of a double chromodomains (Chromatin Organization Modifier) and SNF2-related helicase/ATPase domains. The CHD1/CHD2 sub-family of CHD proteins also contain a C-terminal DNA binding domain that recognizes AT-rich sequences. CHD proteins alter gene expression possibly by modification of chromatin structure thus altering access of the transcriptional apparatus to its chromosomal DNA template. CHD2 is required for normal development of mouse embryos.

Immunogen: This CHD2 antibody (pAb) was raised against a peptide corresponding to the C-terminus of human CHD2.

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:

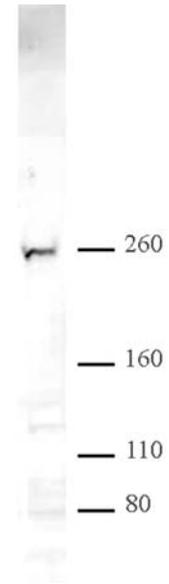
Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 dilution

The addition of 0.1% Tween 20 in all blocking solutions may 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.



CHD2 pAb tested by Western blot.

Detection of CHD2 by Western blot analysis. Nuclear extract of HEK293 cells (25 µg) was probed with CHD2 pAb at a 1:250 dilution.