

## CUL4B antibody (pAb)

**Catalog Nos:** 61445, 61446

**RRID:** AB\_2793640

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 110 kDa

**Background:** **CUL4B** (Cullin 4B) is a core component of multiple cullin-RING-based E3 ubiquitin-protein ligase complexes which mediate the ubiquitination and subsequent proteasomal degradation of target proteins. The functional specificity of the E3 ubiquitin-protein ligase complex depends on the variable substrate recognition subunit. CUL4B may act within the complex as a scaffold protein, contributing to catalysis through positioning of the substrate and the ubiquitin-conjugating enzyme. Plays a role as part of the E3 ubiquitin-protein ligase complex in polyubiquitination of CDT1, Histone H2A, Histone H3 and Histone H4 in response to radiation-induced DNA damage and may act as a transcriptional corepressor by ubiquitination of Histone H2AK119 and recruiting PRC2 complexes.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human CUL4B.

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

### Application Notes:

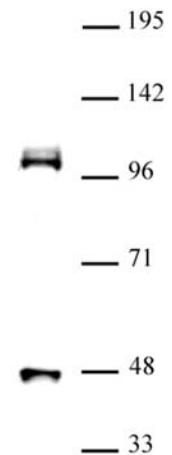
Applications Validated by Active Motif:

WB\*: 1:500 - 1:1,000 dilution

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

**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.



### CUL4B (pAb) tested by Western blot.

Detection of CUL4B by Western blot analysis. Nuclear extract of MCF-7 cells (20 µg) was probed with CUL4B (pAb) at a 1:500 dilution.