

## CTBP2 antibody (pAb)

**Catalog Nos:** 61261, 61262

**RRID:** AB\_2793573

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq, WB

**Reactivity:** Human, Mouse

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 50 kDa

**Background:** The CTBP2 (C-terminal binding protein 2) gene encodes a protein that is a corepressor targeting diverse transcription regulators. It functions in brown adipose tissue (BAT) differentiation. Isoform 2 probably acts as a scaffold for specialized synapses. Isoform 2 is also phosphorylated upon DNA damage, probably by ATM or ATR at Thr179; Ser181 and Ser185. Phosphorylation by HIPK2 on Ser428 induces proteasomal degradation.

**Immunogen:** This CTBP2 antibody was raised against a peptide in the C-terminal region of human CTBP2. This antibody recognizes both isoform 1 and 2.

**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:

ChIP-Seq: 4  $\mu$ g each

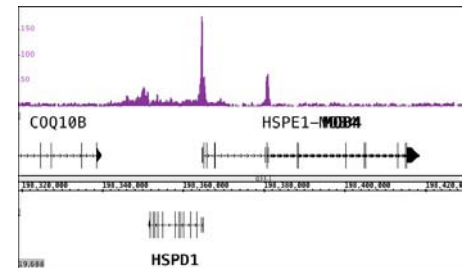
ChIP: 10  $\mu$ l per ChIP

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.

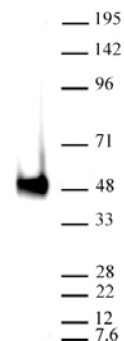
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not 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.



### CTBP2 antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30  $\mu$ g of chromatin from mouse liver and 10  $\mu$ l of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 15 million sequence tags were mapped to identify CTBP2 binding sites. The image shows binding across a regions of chromosome 11. You can view the complete data set in the UCSC Genome Browser



### CTBP2 antibody (pAb) tested by Western blot

Nuclear extract of HeLa cells (10  $\mu$ g) probed with CTBP2 antibody (1:1000).