

CTBP antibody (pAb)

Catalog Nos: 39847, 39848

RRID: AB_2793366

Isotype: Serum

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 50 kDa

Background: CTBP (C-terminal Binding Protein, CTBP1) is a transcriptional repressor that was originally identified by virtue of its ability to bind the C-terminus of E1A proteins. CTBP serves as a co-repressor for a number of DNA binding transcription factors. CTBP is part of a variety of co-repressor complexes, many of which include HDAC proteins. CTBP also interacts with vertebrate Polycomb proteins including Pc2, suggesting that repression of many CTBP-regulated genes also involves recruitment of the Polycomb complex.

Immunogen: This CTBP antibody was raised against a peptide within the C-terminal region of human CTBP.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

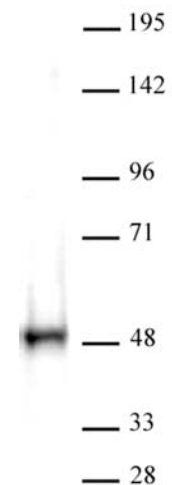
Applications Validated by Active Motif:

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

*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.



CTBP antibody (pAb) tested by Western blot.
Nuclear extract of Jurkat cells (30 µg per lane) probed with CTBP antibody (pAb) at a dilution of 1:2,000.