

CtIP antibody (mAb)

Catalog Nos: 61141, 61942, 61142

RRID: AB_2714164

Clone: 14-1

Isotype: IgG1

Application(s): ChIP, ICC, IF, IP, WB

Reactivity: Human

Quantities: 100 µg, 50 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 120 kDa

Background: CtIP (CTBP interacting protein, RBBP8, RIM, COM1) is a transcriptional regulator and tumor suppressor first identified through its interaction with CTBP (C-Terminal Binding Protein). CtIP binds to the Rb protein and serves as a transcriptional repressor, through recruitment of CtBP and also by interaction with general transcription factors TFIID and TFIIB. CtIP also interacts with Rb-related protein p130 and Ikaros, a master regulator in lymphocyte development. CtIP is phosphorylated upon DNA damage (likely by ATM and/or ATR), interacts with BRCA1 and plays a critical role in facilitating DNA damage repair. Mutations that abolish the interaction between CtIP and BRCA1 result in loss of cell cycle control leading to the development of cancer.

Immunogen: This CtIP antibody was raised against a recombinant protein corresponding to amino acid residues 620-897 of human CtIP.

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

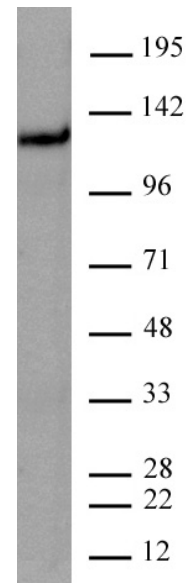
Application Notes:

Applications Validated by Active Motif:

WB: 1 - 2 µg/ml

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.



CtIP antibody (mAb) tested by Western blot
20 µg of HEK293 whole cell extract was run on SDS-PAGE and probed with antibody at 1 µg/ml.