
CNBP Polyclonal Antibody

(Catalog # A70465)

Background

This gene encodes a nucleic-acid binding protein with seven zinc-finger domains. The protein has a preference for binding single stranded DNA and RNA. The protein functions in cap-independent translation of ornithine decarboxylase mRNA, and may also function in sterol-mediated transcriptional regulation. A CCTG expansion from <30 repeats to 75-11000 repeats in the first intron of this gene results in myotonic dystrophy type 2. Multiple transcript variants encoding different isoforms have been found for this gene.

Description

CNBP Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P62633

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 100 to the C-terminus of human CNBP (NP_001120668.1).

Storage

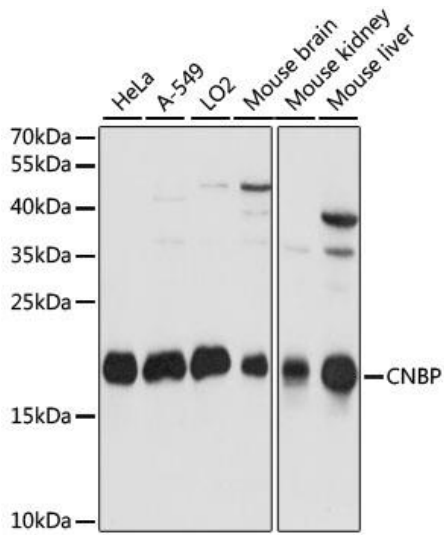
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

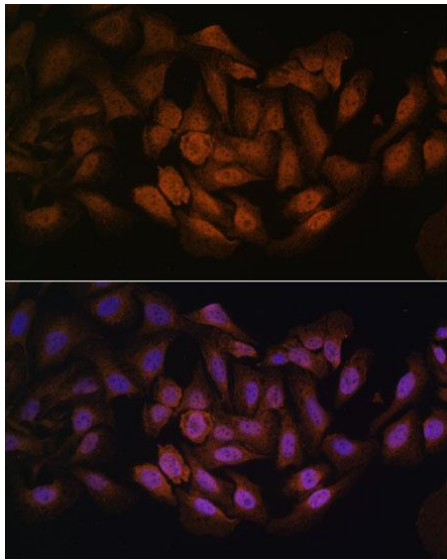
CNBP; CNBP1; DM2; PROMM; RNF163; ZCCHC22; ZNF9; cellular nucleic acid-binding protein

Application

WB, IF; Recommended dilution: WB: 1:500-1:2000, IF:1:50-1:200



Western blot analysis of extracts of various cell lines, using CNBP Polyclonal Antibody at 1:1000 dilution.
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 5s.



Immunofluorescence analysis of U-2 OS cells using CNBP Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.