

ZEB1 antibody (pAb)

Catalog Nos: 61119, 61120

RRID: AB_2793514

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 200 kDa

Background: ZEB1 (Zinc finger E-box-binding homeobox 1, NIL-2-A zinc finger protein Negative regulator of IL2, NIL-2-A, Transcription factor 8, TCF-8) is a POU domain DNA-binding transcriptional regulator that inhibits the expression of E-cadherin and other epithelial genes central to adhesion and epithelial polarity. ZEB1 induces an epithelial-mesenchymal transition (EMT) by recruiting SMARCA4/BRG1. ZEB1 positively regulates neuronal differentiation and it represses RCOR1 transcription activation during neurogenesis. Defects in ZEB1 are the cause of posterior polymorphous corneal dystrophy type 3 (PPCD3).

Immunogen: This ZEB1 antibody was raised against a peptide derived from the N-terminal region of human ZEB1.

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:

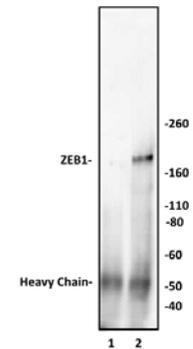
IP: 10 µl per IP

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.

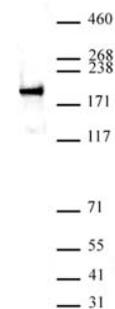
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



ZEB1 antibody (pAb) tested by Immunoprecipitation.

10 µl of ZEB1 antibody was used to immunoprecipitate ZEB1 from 250 µg of HeLa nuclear cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the ZEB1 antibody at a dilution of 1:1,000.



ZEB1 antibody (pAb) tested by Western blot.

Nuclear extract of HeLa cells (40 µg) probed with ZEB1 antibody (1:1,000).