

SIN3A antibody (pAb)

Catalog Nos: 39865, 39866

RRID: AB_2793373

Isotype: Serum

Application(s): ChIP, WB

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 155 kDa

Background: SIN3A (Switch independent 3 homologue) is the mammalian homologue of the yeast protein Sin3. SIN3A is transcriptional co-repressor that is recruited to promoters by sequence-specific DNA binding proteins, such as REST, MXI1 and MAD-MAX. It is part of a large multi-subunit repressor protein complex that also contains HDAC1, HDAC2 and SMRT/NCoR1.

Immunogen: This SIN3A antibody was raised against a peptide derived from the N-terminus of human SIN3A.

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

Application Notes:

Applications Validated by Active Motif:

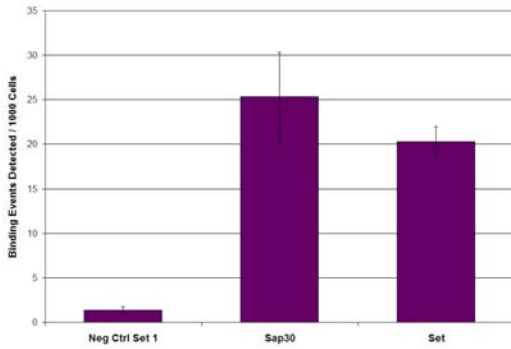
ChIP: 5 - 10 µl per ChIP

WB*: 1:1,000 - 1:5,000 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.

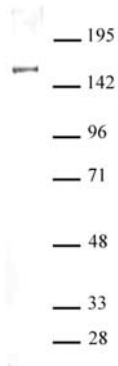
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



SIN3A antibody (pAb) tested by ChIP.

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from TM4 Sertoli cells and 5 µl of SIN3A antibody. ChIP DNA was used in qPCR with the control primer pairs or gene-specific primer pairs as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.



SIN3A antibody (pAb) tested by Western blot.

Nuclear extract of HeLa cells (20 µg) probed with SIN3A antibody at a dilution of 1:5,000.