

## DMAP1 antibody (pAb)

**Catalog Nos:** 61675, 61676

**RRID:** AB\_2793730

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 60 kDa

**Background:** DMAP1 (DNA Methyltransferase 1 Associated Protein 1) is involved in transcription repression and activation. Its interaction with HDAC2 may provide a mechanism for histone deacetylation in heterochromatin following replication of DNA at late firing origins. Can also repress transcription independently of histone deacetylase activity. May specifically potentiate DAXX-mediated repression of glucocorticoid receptor-dependent transcription. Component of the NuA4 histone acetyltransferase (HAT) complex which is involved in transcriptional activation of select genes principally by acetylation of nucleosomal histones H4 and H2A. This modification may both alter nucleosome - DNA interactions and promote interaction of the modified histones with other proteins which positively regulate transcription. This complex may be required for the activation of transcriptional programs associated with oncogene and proto-oncogene mediated growth induction, tumor suppressor mediated growth arrest and replicative senescence, apoptosis, and DNA repair. NuA4 may also play a direct role in DNA repair when recruited to sites of DNA damage. Participates in the nuclear localization of URI1 and increases its transcriptional corepressor activity.

**Immunogen:** This antibody was raised against a peptide within the C-terminal region of human DMAP1.

**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:

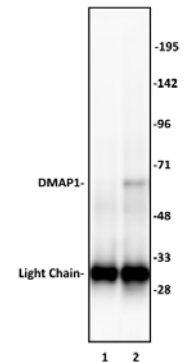
IP: 10 µl per IP

WB\*: 1:500 - 1:2,000 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.

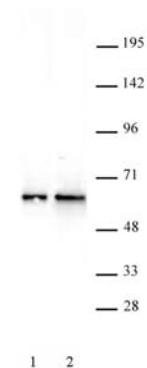
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### DMAP1 antibody (pAb) tested by Immunoprecipitation.

10 µl of DMAP1 antibody was used to immunoprecipitate DMAP1 from 500 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 DMAP1 antibody at a dilution of 1:500.



### DMAP1 antibody (pAb) tested by Western blot.

Nuclear extract (30 µg) of Caco-2 (Lane 1) and HeLa cells (Lane 2) probed with DMAP1 antibody at a 1:500 dilution.