

BAP1 antibody (pAb)

Catalog Nos: 61501, 61502

RRID: AB_2793661

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

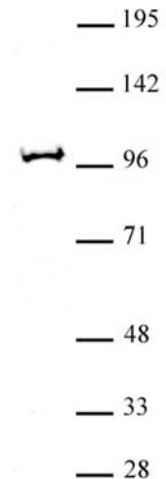
Purification: Affinity Purified

Host: Rabbit

Concentration: 1 μ g/ μ l

Molecular Weight: 100 kDa

Background: **BAP1** (BRCA1 Associated Protein-1 (Ubiquitin Carboxy-Terminal Hydrolase)) is a deubiquitinating enzyme that plays a key role in chromatin by mediating deubiquitination of histone H2A and HCFC1. Catalytic component of the PR-DUB complex, a complex that specifically mediates deubiquitination of histone H2A monoubiquitinated at 'Lys-119' (H2AK119ub1). Does not deubiquitinate monoubiquitinated histone H2B. Acts as a regulator of cell growth by mediating deubiquitination of HCFC1 N-terminal and C-terminal chains, with some specificity toward 'Lys-48'-linked polyubiquitin chains compared to 'Lys-63'-linked polyubiquitin chains. Deubiquitination of HCFC1 does not lead to increase stability of HCFC1. Interferes with the BRCA1 and BARD1 heterodimer activity by inhibiting their ability to mediate ubiquitination and autoubiquitination. It however does not mediate deubiquitination of BRCA1 and BARD1. Acts as a tumor suppressor.



Immunogen: This antibody was raised against a peptide within the C-terminal region of human BAP1.

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

BAP1 antibody (pAb) tested by Western blot.
Nuclear extract of P19 cells (25 μ g) probed with the BAP1 antibody (1:1,000 dilution).

Application Notes:

Applications Validated by Active Motif:

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. The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by 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.