

## ATAD2 antibody (pAb)

**Catalog Nos:** 61369, 61370

**RRID:** AB\_2793609

**Isotype:** IgG

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

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** .46  $\mu$ g/ $\mu$ l

**Molecular Weight:** 160 kDa

**Background:** **ATAD2** (ATPase family AAA domain-containing protein 2) may be a transcriptional coactivator of the nuclear receptor ESR1 required to induce the expression of a subset of estradiol target genes, such as CCND1, MYC and E2F1. May play a role in the recruitment or occupancy of CREBBP at some ESR1 target gene promoters. May be required for histone hyperacetylation. Involved in the estrogen-induced cell proliferation and cell cycle progression of breast cancer cells.

**Immunogen:** This antibody was raised against a peptide within the C-terminal region of human ATAD2.

**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: 5 - 10  $\mu$ l per IP

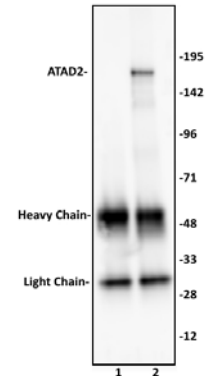
WB\*: 1:500 - 1:2,000 dilution

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. Individual optimization may be required.

\*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.

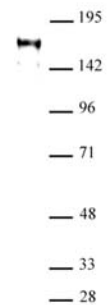
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not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



### ATAD2 antibody (pAb) tested by Immunoprecipitation.

10  $\mu$ l of ATAD2 antibody was used to immunoprecipitate ATAD2 from 250  $\mu$ g of Jurkat nuclear cell extract (lane 2). 10  $\mu$ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the ATAD2 antibody at a dilution of 1:1,000.



### ATAD2 antibody (pAb) tested by Western blot.

ATAD2 detection by Western blot. The analysis was performed using 40  $\mu$ g Jurkat nuclear cell extract and CENP-B at a 1:1,000 dilution.