

## ATRX antibody (pAb)

**Catalog Nos:** 61049, 61050

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 0.85 µg/µl

**Molecular Weight:** 300 kDa

**Background:** ATRX (alpha thalassemia/mental retardation syndrome X-linked protein, X-linked helicase II, XNP) is a transcriptional regulator and a member of the SWI/SNF family of chromatin remodeling proteins. It contains an SNF2 domain and helicase domain, and a PHD finger. ATRX undergoes cell cycle-dependent phosphorylation, which regulates its nuclear matrix and chromatin association, and suggests its involvement in interphase gene regulation and chromosomal segregation in mitosis. Mutations of the ATRX gene are associated with an X-linked mental retardation syndrome most often accompanied by alpha-thalassemia (ATRX) syndrome. These mutations have been shown to cause diverse changes in the pattern of DNA methylation, which may provide a link between chromatin remodeling, DNA methylation, and gene expression in developmental processes.

**Immunogen:** This ATRX antibody was raised against a peptide within the C-terminus of human ATRX.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

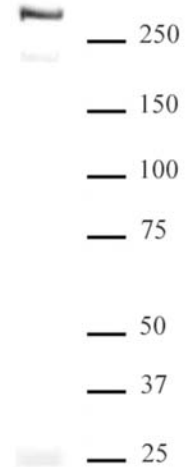
### Application Notes:

Applications Validated by Active Motif:

WB: 1:1,000 - 1:5,000 dilution

**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.



### Western blot of ATRX antibody.

Nuclear extract of CCF-STGG1 cells (40 µg per lane) probed with ATRX antibody (pAb) at a dilution of 1:2,500.