

## TRIM37 antibody (pAb)

**Catalog Nos:** 61747, 61748

**RRID:** AB\_2793755

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human, Mouse

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 120 kDa

**Background:** **TRIM37** (Tripartite Motif Containing 37) is an E3 ubiquitin-protein ligase required to prevent centriole reduplication (PubMed:15885686, PubMed:23769972). Probably acts by ubiquitinating positive regulators of centriole reduplication (PubMed:23769972). Mediates monoubiquitination of Lys-119 of histone H2A (H2AK119Ub), a specific tag for epigenetic transcriptional repression: associates with some Polycomb group (PcG) multiprotein PRC2-like complex and mediates repression of target genes (PubMed:25470042). Has anti-HIV activity (PubMed:24317724).

**Immunogen:** This antibody was raised against a peptide within the C-terminal region of human TRIM37.

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

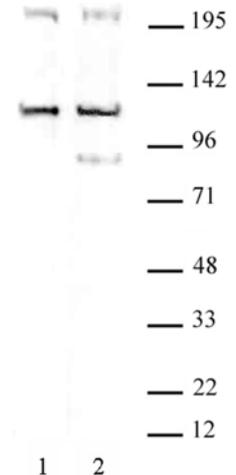
### Application Notes:

Applications Validated by Active Motif:

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



### TRIM37 antibody (pAb) tested by Western blot.

TRIM37 antibody detection by Western blot. The analysis was performed using 30 µg of human MDA-MB-231 nuclear cell extract (Lane 1) or mouse F9 nuclear extract (Lane 2) and TRIM37 antibody at a 1:500 dilution.