

## TRIM28 antibody (mAb)

**Catalog Nos:** 61173, 61174

**RRID:** AB\_2793537

**Clone:** 1Tb-1A9

**Isotype:** IgG2b

**Application(s):** ChIP, ICC, IF, WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 90 kDa

**Background:** Tripartite Motif-containing Protein 28 (TRIM28), also known as KAP1 (KRAB Associated Protein 1, and TIF1β, Transcriptional Intermediary factor 1β, is an important co-repressor protein that plays a critical role in proliferation and differentiation of both normal and tumor cells. It functions as a co-repressor for KRAB domain-containing zinc finger proteins (KRAB-ZFPs) and mediates gene silencing by recruiting CHD3, a subunit of the Nucleosome Remodeling and Deacetylation (NuRD) complex. Part of its mechanism involves coordinating the increase in H3K9 methylation at promoters of silenced genes, potentially through direct recruitment of SETDB1, an H3K9 methyltransferase involved in the formation of heterochromatin. TRIM28 is also involved in maintaining pluripotency and is required for terminal differentiation of embryonic stem cells. TRIM28 protein levels are increased in liver, gastric, lung, breast, and prostate cancer, and gastric cancer patients with high levels of TRIM28 show a significantly lower survival rate. TRIM28 contains two zinc fingers and a PHD finger.

**Immunogen:** This TRIM28 antibody was raised against a full-length protein corresponding to mouse TRIM28.

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

### Application Notes:

Validated Applications:

ICC/IF: 0.5 µg/ml dilution

Published Applications:

ChIP

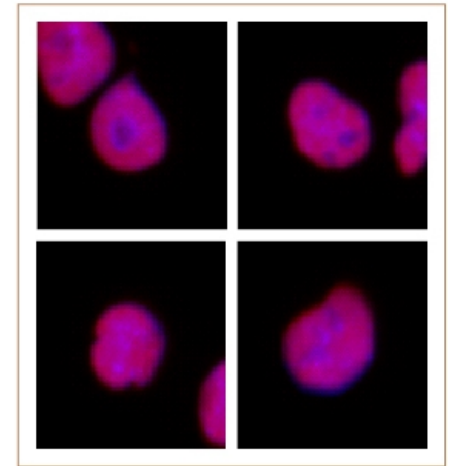
ICC/IF

WB

See references for more information. Individual optimization may be required.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### TRIM28 antibody (mAb) tested by Immunofluorescence.

Formaldehyde fixed HeLa cells stained with TRIM28 antibody at a 0.5 µg/ml dilution.