

## Tada2a antibody (mAb)

---

**Catalog Nos:** 61333, 61334

**RRID:** AB\_2614995

**Clone:** 2AD-2A1

**Isotype:** IgG1k

**Application(s):** IP

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 0.77 µg/µl

**Molecular Weight:** 54 kDa

---

**Background:** **Tada2a** (transcriptional adaptor 2A) is a component of the ATAC complex, a complex with histone acetyltransferase activity on histones H3 and H4. Required for the function of some acidic activation domains, which activate transcription from a distant site (By similarity). Binds double-stranded DNA. Binds dinucleosomes, probably at the linker region between neighboring nucleosomes. Plays a role in chromatin remodeling.

**Immunogen:** This antibody was raised against a synthetic peptide corresponding to amino acid residue(s) 1-17 of human Tada2a.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide.

**Application Notes:**

Published Applications:

IP

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.