

Trrap antibody (mAb)

Catalog No: 61335

RRID: AB_2793596

Clone: 2TRR-1B3

Isotype: IgG2a

Application(s): ChIP, IP

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 0.47 µg/µl

Molecular Weight: 450 kDa

Background: Trrap (transformation/transcription domain-associated protein) is an adapter protein, which is found in various multiprotein chromatin complexes with histone acetyltransferase activity (HAT), which gives a specific tag for epigenetic transcription activation. Component of the NuA4 histone acetyltransferase complex which is responsible for acetylation of nucleosomal histones H4 and H2A. Plays a central role in MYC transcription activation, and also participates in cell transformation by MYC. Required for p53/TP53-, E2F1- and E2F4-mediated transcription activation. Also involved in transcription activation mediated by the adenovirus E1A, a viral oncoprotein that deregulates transcription of key genes. Probably acts by linking transcription factors such as E1A, MYC or E2F1 to HAT complexes such as STAGA thereby allowing transcription activation. Probably not required in the steps following histone acetylation in processes of transcription activation. May be required for the mitotic checkpoint and normal cell cycle progression.

Immunogen: This antibody was raised against a synthetic peptide corresponding to amino acid residue(s) 2023-2042 of human Trrap.

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:

ChIP

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.