

NAT10 antibody (pAb)

Catalog Nos: 65703, 65903, 65704**RRID:** AB_3216304**Application(s):** WB**Reactivity:** Human**Quantities:** 100 µg, 50 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 130 kDa

Background: NAT10 (N-Acetyltransferase 10) antibody is a RNA cytidine acetyltransferase that catalyzes the formation of N(4)- acetylcytidine (ac4C) modification on mRNAs, 18S rRNA and tRNAs and mediates the formation of ac4C at position 1842 in 18S rRNA. NAT10 requires the tRNA-binding adapter protein THUMP1 for full tRNA acetyltransferase activity but not for 18S rRNA acetylation. In addition to RNA acetyltransferase activity, NAT10 is also able to acetylate lysine residues of proteins, such as histones, microtubules, p53/TP53 and MDM2, in vitro. It activates telomerase activity by stimulating the transcription of TERT, and may also regulate telomerase function by affecting the balance of telomerase subunit assembly, disassembly, and localization. NAT10 involved in the regulation of centrosome duplication by acetylating CENATAC during mitosis, promoting SASS6 proteasome degradation.

Immunogen: This antibody was raised against a human RNA cytidine acetyltransferase isoform a.

Buffer: Purified IgG supplied in 62 mM Trisodium citrate, 87 mM Tris pH 7, 30% glycerol, 0.035 % Sodium Azide.

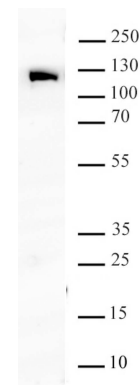
Application Notes:

Applications Validated by Active Motif:

WB: 1 - 4 µg 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.

**NAT10 antibody tested by Western blot.**

K562 Nuclear Extract (20 µg per lane) was probed with NAT10 antibody (1 µg/mL dilution).