

INHAT-1/TAF-1 α /TAF-1 β antibody (pAb)**Catalog No:** 39202**RRID:** AB_2793183**Application(s):** ICC, WB**Reactivity:** Human, Mouse, Rat**Volume:** 200 μ l**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Concentration:** 0.5 μ g/ μ l**Molecular Weight:** 37 kDa

Background: INHAT (also designated TAF-1 α , TAF-1 β and Inhibitor of Histone Acetyltransferases complex) binds to histones and masks them from being substrates for histone acetyltransferases (HATs). Endogenous INHAT subunits, which include the Set/TAF-1 β oncoprotein, and pp32 associate with chromatin *in vivo* and can block coactivator-mediated transcription when transfected in cells. Single phosphorylations within the Histone H3 tail blocks binding of INHAT, as well as simultaneous acetylation of multiple lysine residues.

Immunogen: This INHAT-1/TAF-1 α /TAF-1 β antibody was raised against a mixture of synthetic peptides corresponding to amino acid residues 66-81 and 135-151 of human INHAT-1/TAF-1 α /TAF-1 β .

Buffer: PBS containing 0.02% sodium azide and 0.1 mg/ml BSA. Sodium azide is highly toxic.

Application Notes:

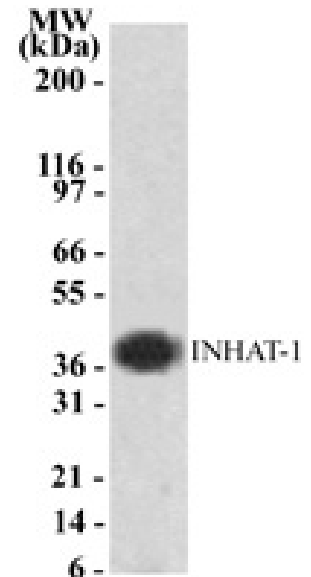
Applications Validated by Active Motif:

WB: 1 - 2 μ g/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. 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.

**INHAT-1/TAF-1 α /TAF-1 β pAb tested by Western blot.**

INHAT-1 detection by Western blot. The analysis was performed using Jurkat whole-cell extract and INHAT-1/TAF-1 α /TAF-1 β pAb at a 2 μ g/ml dilution.