

Supt3h antibody (mAb)

Catalog Nos: 61343, 61344

RRID: AB_2614996

Clone: 1SP-2C10

Isotype: IgG1k

Application(s): ChIP, IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 38 kDa

Background: Supt3h (suppressor of Ty 3 homolog (*S. cerevisiae*)) is a probable transcriptional activator and functions as a component in multiple histone acetyltransferase (HAT) complexes such as STAGA (SPT3-TAF-GCN5 acetylase), PCAF (p300/CBP associated factor), and TFTC (TBP-free TAF complex).

Immunogen: This antibody was raised against a peptide corresponding to amino acid residue(s) 10-27 of human Supt3h.

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

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.