

---

## Histone Acetyltransferase KAT5 Polyclonal Antibody

(Catalog #A54421)

---

### Background

Catalytic subunit of the NuA4 histone acetyltransferase complex which is involved in transcriptional activation of select genes principally by acetylation of nucleosomal histones H4 and H2A. This modification may both alter nucleosome-DNA interactions and promote interaction of the modified histones with other proteins which positively regulate transcription. This complex may be required for the activation of transcriptional programs associated with oncogene and proto-oncogene mediated growth induction, tumor suppressor mediated growth arrest and replicative senescence, apoptosis, and DNA repair. NuA4 may also play a direct role in DNA repair when recruited to sites of DNA damage. Directly acetylates and activates ATM. In case of HIV-1 infection, interaction with the viral Tat protein leads to KAT5 polyubiquitination and targets it to degradation.

### Description

Histone Acetyltransferase KAT5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q92993

### Purification

Protein G purified

### Immunogen

Recombinant Human Histone acetyltransferase KAT5 protein (3-512AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

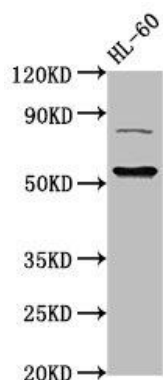
### Alternative Names

60 kDa Tat-interactive protein, Histone acetyltransferase HTATIP, Lysine acetyltransferase 5, cPLA(2)-interacting protein, KAT5, HTATIP, TIP60

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000, IF:1:50-1:500

---



#### Western Blot

Positive WB detected in: HL60 whole cell lysate

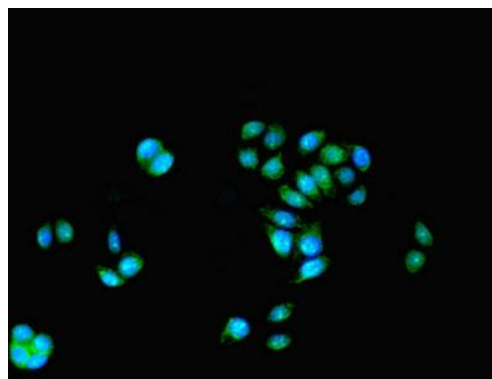
All lanes: KAT5 antibody at 3µg/ml

Secondary

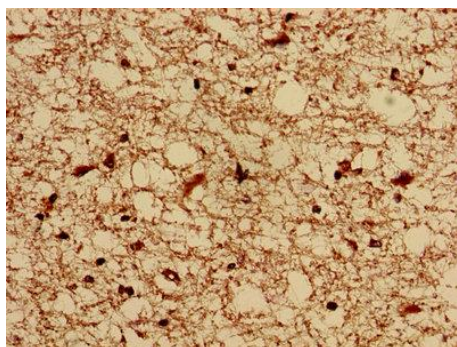
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 59, 54, 62, 57 kDa

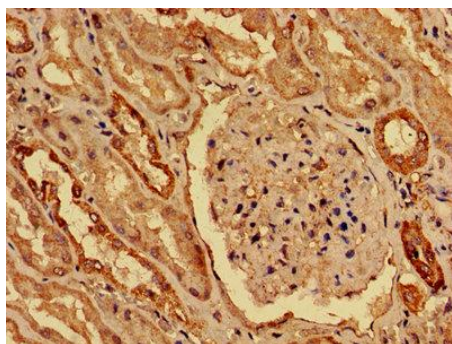
Observed band size: 59 kDa



Immunofluorescent analysis of PC-3 cells using KAT5 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



IHC image of KAT5 Antibody diluted at 1:800 and staining in paraffin-embedded human brain tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of KAT5 Antibody diluted at 1:800 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.