

# **GCN5** Polyclonal Antibody

(Catalog #A-4013)

## **Background**

GCN5, a putative transcriptional adapter in humans and yeast, possesses histone acetyltransferase (HAT) activity. GCN5 HATs operate in multi-subunit complexes (e.g., SAGA and ADA) to acetylate lysine residues in histone tails, preferentially those of histone H3. Also, GCN5 is required for the activation of a number of genes involved in amino acid biosynthesis, respiration, and other pathways controlled by the SWI/SNF gene products.

## **Description**

GCN5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

#### **Formulation**

Liquid. PBS containing 50% glycerol, 0.5% BSA\* and 0.02% sodium azide.

## Specificity

Human

# Isotype

IgG

#### **Uniprot ID**

Q92830

#### **Purification**

Affinity Purified

## **Immunogen**

Synthesized peptide derived from the C-terminal region of human GCN5.

#### Storage

Store at -20°C or -80°C. Avoid repeated freeze.

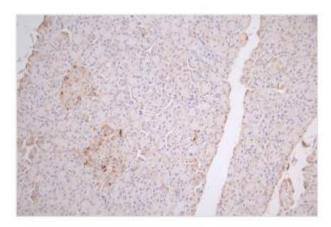
#### **Alternative Names**

KAT2A; GCN5; GCN5L2; HGCN5; Histone acetyltransferase KAT2A; General control of amino acid synthesis protein 5-like 2; Histone acetyltransferaseGCN5; HsGCN5; Lysine acetyltransferase 2A; STAF97

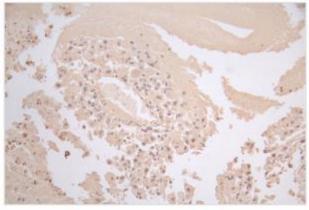
# **Application**

IF, IHC, ELISA; Recommended Dilution; IF:1:20-1:300, IHC:1:100-1:300, ELISA:1:20000

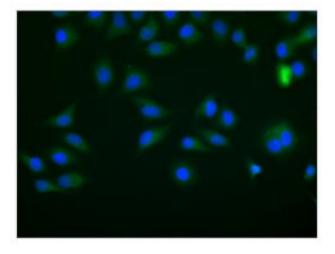
\*The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.



IHC image of GCN5 Polyclonal Antibody diluted at 1:66 and staining in paraffin-embedded human pancreatic 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 Goat anti-rabbit polymer IgG labeled by HRP and visualized using 0.05% DAB.



IHC image of GCN5 Polyclonal Antibody diluted at 1:66 and staining in paraffin-embedded human melanoma cancer 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 Goat anti-rabbit polymer IgG labeled by HRP and visualized using 0.05% DAB.



Immunofluorescence staining of Hela cell with GCN5 Polyclonal Antibody at 1:30, counter-stained with DAPI. The cells were fixed in 4% formaldehyde and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4C. The secondary antibody was Alexa Fluor 488-congugated AffiniPure Goat Anti-Rabbit IgG(H+L).