

## 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 multisubunit 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, Mouse

### Isotype

IgG

### Uniprot ID

Q92830

### Purification

Affinity Purified

### Immunogen

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

### Storage

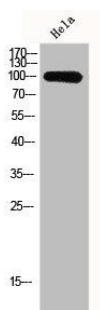
Store at -20°C or -80°C. Avoid repeated freeze.

### Alternative Names

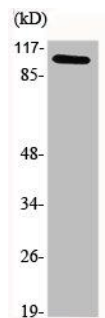
KAT2A; GCN5; GCN5L2; HGCN5; Histone acetyltransferase KAT2A; General control of amino acid synthesis protein 5-like 2; Histone acetyltransferase GCN5; HsGCN5; Lysine acetyltransferase 2A; STAF97

### Application

WB, IHC, ELISA; Recommended dilution: WB:1/500-1/2000, IHC:1/100-1/300, ELISA:1/20000



Western Blot analysis of HeLa cells using GCN5 Polyclonal Antibody



Western Blot analysis of various cells using GCN5 Polyclonal Antibody