

GCN5 antibody (mAb)

Catalog No: 39975

RRID: AB_2793415

Clone: 2GC-2C11

Isotype: IgG1

Application(s): ICC, IF, WB

Reactivity: Human

Quantity: 100 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 0.48 µg/µl

Molecular Weight: 95 kDa

Background: GCN5 (General control of nutrition 5, KAT2A) is a histone acetyltransferase (HAT) and serves as a transcriptional activator. GCN5 is a component of the SAGA and ATAC complexes, which have histone acetyltransferase activity for histones H3 and H4. GCN5 is required for neural tube closure and embryonic development.

Immunogen: This GCN5 antibody was raised against a synthetic peptide corresponding to amino acids 118-132 of human GCN5.

Buffer: PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Published Applications:

ICC/IF

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.