

GCN5 Bromodomain

Catalog No. 15-0026
Lot No. 13228001
Pack Size 100 µg

Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 36 kDa **Epitope Tag** GST

Product Description:

Recombinant human GCN5 Bromodomain (KAT2A, GCN5L2, HGCN5, accession number Q92830, amino acids 721 to 837) containing an N-terminal GST tag, expressed in *E. coli*. GCN5 is an histone acetyltransferase that catalyzes the acetylation of multiple sites on H3 and H4. GCN5 is a component of several protein complexes involved in transcriptional activation including SAGA and ATAC. The bromodomain for GCN5 recognizes multiple sites of acetylation on all four core histones.

Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

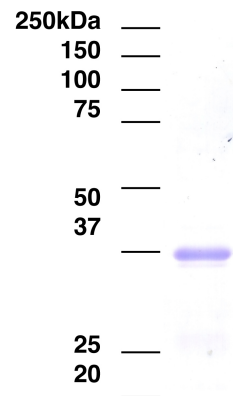
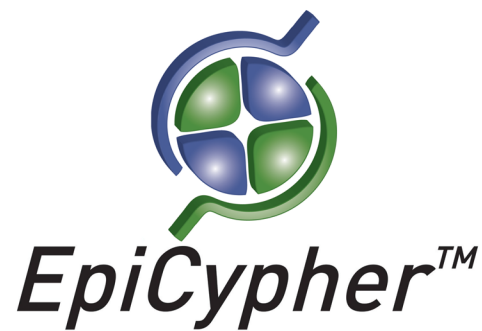
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Recombinant GCN5 Bromodomain is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: GCN5 Bromodomain (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.