
GCN5 Protein

(Catalog No. E24001)

Description

Human GCN5, also known as KAT2A, GCNL2, or PCAF- β , GenBank Accession No. NM_021078, a.a. 362-837(end), with C- terminal FLAG-tag.

Molecular Weight (kDa)

55 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 μ g/ml FLAG peptide, and 20% glycerol.

Source

Expressed in E. coli

Size

100 μ g

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Glutathione Sepharose

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
GCN5 Protein	100 μ g	E24001-1

This product is for research purposes only. Not intended for use in diagnostic procedures.