

## Recombinant KAT7 protein

---

**Catalog No:** 31489, 31889

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.6 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant KAT7 protein is supplied at a concentration of 0.6 µg/µl in 25 mM Hepes pH 7.5, 300 mM NaCl, 5% Glycerol, 0.04% Triton X-100, 0.2 mM TCEP, 0.2 mg/ml 3× Flag peptide.

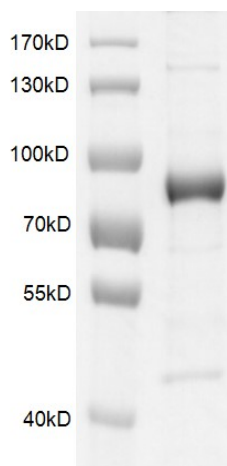
**Background:** **KAT7** (K(Lysine) Acetyltransferase 7) is a MYST-family histone acetyltransferase (HAT) that forms a complex with ING4/5, JADE1/2/3, and Eaf6, which is found enriched on coding regions of genes, supporting a role in transcriptional elongation. In addition, KAT7 is thought to participate in events controlling the initiation of DNA replication.

**Protein Details:** Full length recombinant KAT7 protein (accession number NP\_001186084.1) was expressed in Sf9 cells and contains an N-terminal Flag tag with a molecular weight of 67.2 kDa. The recombinant protein is >85% pure by SDS-PAGE.

**Application Notes:** Recombinant KAT7 protein is suitable for use in enzyme kinetics, inhibitor screening, and selectivity profiling.

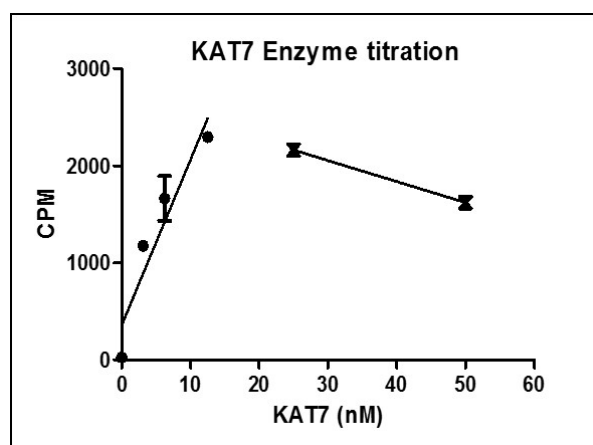
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



#### Recombinant KAT7 protein gel.

Recombinant KAT7 was run on an 8% SDS-PAGE gel and stained with Coomassie blue.



#### Recombinant KAT7 protein activity assay

Recombinant KAT7 activity titration was assessed using a Radiometry assay. Titration curves were generated to show signal response in the presence of histone peptide substrate at increasing KAT7 concentrations. This data was generated and kindly provided courtesy of ChemPartner.