

Recombinant KAT7 protein

Catalog No: 31489, 31889

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.4 µg/µl

Source: Human

Buffer Contents: Full length recombinant KAT7 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

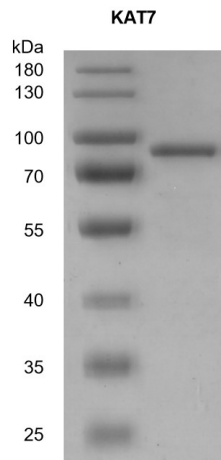
Background: KAT7 (K(Lysine) Acetyltransferase 7) is a MYST-family histone acetyltransferase (HAT) that forms a complex with ING4/5, JADE1/2/3, and Eaf6, which is found enriched on coding regions of genes, supporting a role in transcriptional elongation. In addition, KAT7 is thought to participate in 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 >90% pure by SDS-PAGE.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

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

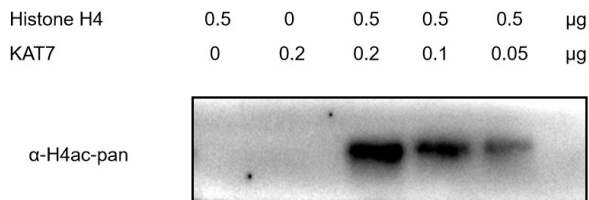
10% SDS-PAGE Coomassie staining

MW: 67.2 kDa

Purity: $\geq 95\%$

Recombinant KAT7 protein activity assay

0.5 μg of Histone H4 (Cat# 31493) was incubated with 0 μg , 0.05 μg , 0.1 μg , and 0.2 μg KAT7 protein in 20 μl reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X-100, 2 mM MgCl_2 , 1 mM TCEP, 20 μM acetyl-CoA for 2 hr at room temperature. 2.5 μl of reaction products were run on a 12.5% SDS-PAGE and detected by H4ac-pan antibody (Cat# 39243).



Western Blot was used for activity detection.