

## Recombinant KAT6B / MORF (718-1008) protein

---

**Catalog No:** 81224, 81924

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

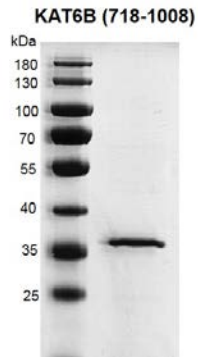
**Buffer Contents:** Recombinant KAT6B / MORF (718-1008) protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

**Background:** KAT6B (Lysine Acetyltransferase 6B), also called as MOZ2, MORF or MYST4, is a histone acetyltransferase and component of the MOZ/MORF protein complex. It has a histone H3 acetyltransferase activity. In addition to its acetyltransferase activity, the encoded full length protein has transcriptional activation activity in its N-terminal end and transcriptional repression activity in its C-terminal end. So it may be involved in both positive and negative regulation of transcription. KAT6B is necessary for RUNX2-dependent transcriptional activation and could be involved in brain development. Mutations have been found in patients with genitopatellar syndrome. A translocation of KAT6B gene and the CREBBP gene results in acute myeloid leukemias. It may also be involved in cerebral cortex development.

**Protein Details:** Recombinant KAT6B / MORF (718-1008) protein that includes amino acids 718-1008 of human KAT6B protein (accession number NP\_036462.2) was expressed in a baculovirus expression system and contains an N-terminal FLAG tag. The molecular weight of the protein is 35.7 kDa. This product is the catalytic domain of KAT6B protein.

**Application Notes:** Recombinant KAT6B / MORF (718-1008) 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

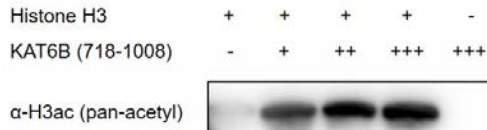


**Recombinant KAT6B / MORF (718-1008) protein gel**

10% SDS-PAGE gel with Coomassie blue staining

MW: 35.7 kDa

Purity: > 93%



**Recombinant KAT6B / MORF (718-1008) protein activity assay**

0.5  $\mu$ g Histone H3.1 (Cat. No. 31296) was incubated with 0  $\mu$ g (-), 0.1  $\mu$ g (+), 0.2  $\mu$ g (++) , 0.4  $\mu$ g (+++) KAT6B / MORF (718-1008) protein in a 20  $\mu$ l reaction system containing 50 mM Tris-HCl pH 8.6, 2 mM MgCl<sub>2</sub>, 1 mM TCEP, 0.02% Triton X-100 and 20  $\mu$ M Acetyl-CoA for 2 hr at RT. Reaction products were detected by Western blot using H3ac (pan-acetyl) antibody (Cat. No. 61637).