

Recombinant KAT6A / MOZ (488-778) protein

Catalog No: 81223, 81923

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.4 µg/µl

Source: Human

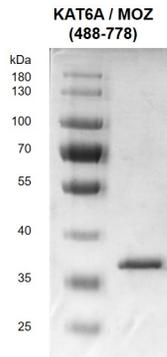
Buffer Contents: Recombinant KAT6A / MOZ (488-778) protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: KAT6A (Lysine Acetyltransferase 6A), also called as MOZ, MYST3 or ZNF220, is a histone acetyltransferase that acetylates lysine residues in histone H3 and histone H4 (in vitro). It is a component of the MOZ / MORF complex which has a histone H3 acetyltransferase activity. KAT6A may act as a transcriptional coactivator for RUNX1 and RUNX2. It also acetylates p53 / TP53 at 'Lys-120' and 'Lys-382' and controls its transcriptional activity via association with PML. This protein is composed of a nuclear localization domain, a double C2H2 zinc finger domain that binds to acetylated histone tails, a histone acetyl-transferase domain, a glutamate/aspartate-rich region, and a serine- and methionine-rich transactivation domain.

Protein Details: Recombinant KAT6A / MOZ (488-778) protein that includes amino acids 488-778 of human KAT6A protein (accession number NP_006757.2) was expressed in a baculovirus expression system and contains an N-terminal FLAG tag. The molecular weight of the protein is 35.5 kDa. This product is the catalytic domain of KAT6A protein.

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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant KAT6A / MOZ (488-778) protein gel

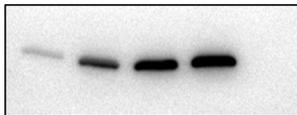
10% SDS-PAGE gel with Coomassie staining

MW: 35.5 kDa

Purity: > 95%

Histone H3.1	+	+	+	+	-
KAT6A	-	+	++	+++	+++

α -H3ac (pan-acetyl)



Recombinant KAT6A / MOZ (488-778) protein activity assay

0.5 μ g Histone H3.1 (Cat. No. 31296) was incubated with 0 μ g (-), 0.1 μ g (+), 0.2 μ g (++) , 0.4 μ g (+++) KAT6A / MOZ (488-778) protein in a 20 μ l reaction system containing 50 mM Tris-HCl pH 8.6, 2 mM MgCl₂, 1 mM TCEP, 0.02% Triton X-100 and 20 μ M Acetyl-CoA for 2 hr at RT. Reaction products were detected by Western blot using H3 pan-acetyl antibody (Cat. No. 61637).