

Recombinant ALKBH4 protein

Catalog No: 81131, 81831

Lot No: 14318001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.3 µg/µl

Source: Human

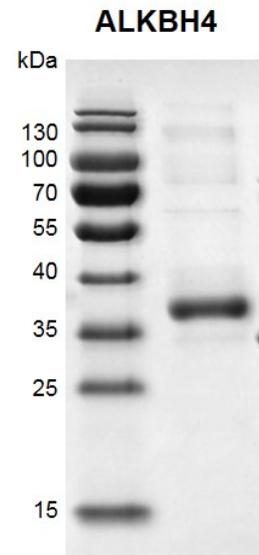
Buffer Contents: Recombinant ALKBH4 protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol, and 0.5 mM TCEP

Background: ALKBH4 (Alpha-Ketoglutarate-Dependent Dioxygenase AlkB Homolog 4) is a member of ALKBH protein family. It's a dioxygenase that mediates demethylation of actin monomethylated at Lys-84 (K84me1), thereby acting as a regulator of actomyosin-processes. Demethylation of actin K84me1 is required for maintaining actomyosin dynamics supporting normal cleavage furrow ingression during cytokinesis and cell migration. ALKBH4 may also be involved in transcription regulation.

Protein Details: Full length ALKBH4 protein (accession number NP_060091.1) was expressed in *E. coli* cells with an N-terminal 6Å~His tag. The molecular weight of ALKBH4 is 37.4 kDa.

Application Notes: Recombinant ALKBH4 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 ALKBH4 protein gel
12.5% SDS-PAGE Coomassie staining
MW: 37.4 kDa
Purity: >85%