

## Recombinant ALKBH1 protein

**Catalog No:** 81128, 81828

**Expressed In:** *E. coli*

**Quantity:** 50, 1000 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

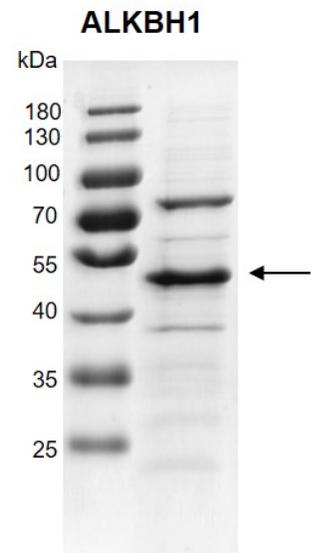
**Buffer Contents:** Recombinant ALKBH1 protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

**Background:** ALKBH1 (AlkB Homolog 2, Alpha-Ketoglutarate Dependent Dioxygenase) is a member of ALKBH protein family. It can repair alkylated DNA and RNA containing 1-methyladenine and 3-methylcytosine by oxidative demethylation. ALKBH1 can also repair alkylated DNA containing 1-ethenoadenine (in vitro). Has strong preference for double-stranded DNA. It can release modified bases from both DNA and RNA. The oxidative demethylation requires molecular oxygen, alpha-ketoglutarate and iron.

**Protein Details:** Full length ALKBH1 protein (accession number NP\_006011.2) was expressed in *E. coli* cells with an N-terminal 6×His-Tag. The molecular weight of ALKBH1 is 47.6 kDa.

**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 this product is shown.

**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 ALKBH1 protein gel**  
10% SDS-PAGE Coomassie staining  
MW: 47.6 kDa  
Purity: >60%