

Recombinant ALDOC protein

Catalog No: 81305, 81605

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 1 µg/µl

Source: Human

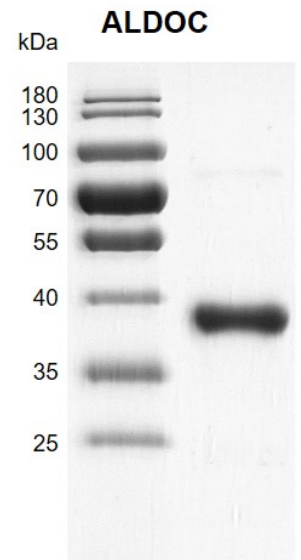
Buffer Contents: Recombinant ALDOC protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: ALDOC (Aldolase, Fructose-Bisphosphate C) is a member of the class I fructose-biphosphate aldolase gene family. Expressed specifically in the hippocampus and Purkinje cells of the brain, the encoded protein is a glycolytic enzyme that catalyzes the reversible aldol cleavage of fructose-1,6-biphosphate and fructose 1-phosphate to dihydroxyacetone phosphate and either glyceraldehyde-3-phosphate or glyceraldehyde, respectively.

Protein Details: Full length human ALDOC protein was expressed in *E. coli* cells (accession number NP_005156.1) with a C-terminal 6×His tag. The molecular weight of the protein is 40.5 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 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 arrival.



Recombinant ALDOC protein gel
12.5% SDS-PAGE with Coomassie blue staining

MW: 40.5 kDa

Purity: >92%