

Recombinant ALDOA protein

Catalog No: 81304, 81604

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 1.8 µg/µl

Source: Human

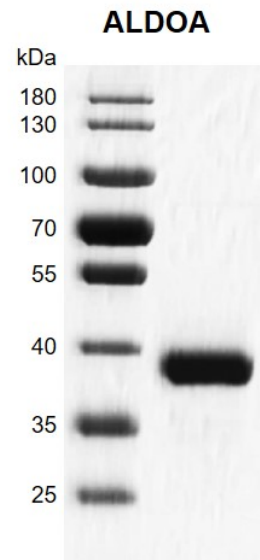
Buffer Contents: Recombinant ALDOA protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

Background: ALDOA (Aldolase, Fructose-Bisphosphate A) is a member of the class I fructose-bisphosphate aldolase protein family. The encoded protein is a glycolytic enzyme that catalyzes the reversible conversion of fructose-1,6- biphosphate to glyceraldehyde 3-phosphate and dihydroxyacetone phosphate. Three aldolase isozymes (A, B, and C), encoded by three different genes, are differentially expressed during development. Mutations in this gene have been associated with Glycogen Storage Disease XII, an autosomal recessive disorder associated with hemolytic anemia. Disruption of this gene also plays a role in the progression of multiple types of cancers.

Protein Details: Full length human ALDOA protein was expressed in *E. coli* cells (accession number NP_908930.1) 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 ALDOA protein gel
10% SDS-PAGE with Coomassie blue staining

MW: 40.5 kDa

Purity: >95%