

Recombinant CTBP1 protein

Catalog No: 81250, 81950

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 3.6 µg/µl

Source: Human

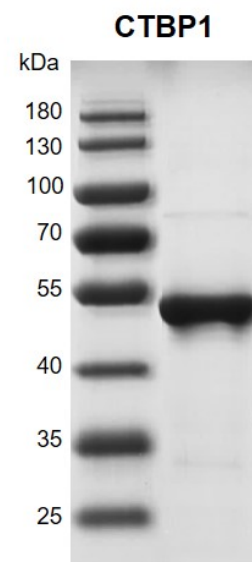
Buffer Contents: Recombinant CTBP1 protein is supplied in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 10% glycerol, 0.5 mM TCEP.

Background: CTBP1 (C-Terminal Binding Protein 1), also written as CtBP1, is a protein that binds to the C-terminus of adenovirus E1A proteins. This phosphoprotein is a transcriptional repressor and may play a role during cellular proliferation. It's a corepressor targeting diverse transcription regulators such as GLIS2 or BCL6. It also has dehydrogenase activity. Involved in controlling the equilibrium between tubular and stacked structures in the Golgi complex. Functions in brown adipose tissue (BAT) differentiation.

Protein Details: Recombinant CTBP1 protein was expressed in *E. coli* as the full length protein (accession number (NP_001319.1) with an N-terminal 6x-His Tag. The molecular weight of the protein is 51.1 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 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 guaranteed for 6 months from date of arrival.



Recombinant CTBP1 protein
10% SDS-PAGE Coomassie staining

MW: 51.1 kDa

Purity: >90%