

Recombinant CHD3 protein

Catalog No: 81386, 81686

Expressed In: Baculovirus

Quantity: 10, 1000 µg

Concentration: 0.25 µg/µl

Source: Human

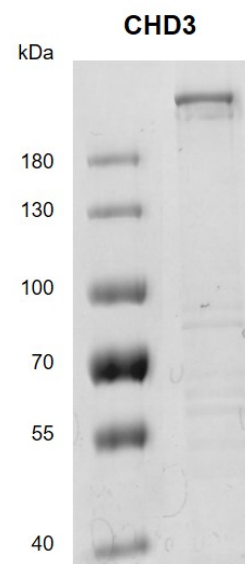
Buffer Contents: Recombinant CHD3 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: CHD3 (Chromodomain Helicase DNA Binding Protein 3) is a member of the CHD family of proteins which are characterized by the presence of chromo (chromatin organization modifier) domains and SNF2-related helicase/ ATPase domains. This protein is one of the components of a histone deacetylase complex referred to as the Mi-2/ NuRD complex which participates in the remodeling of chromatin by deacetylating histones and is required for anchoring centrosomal pericentrin in both interphase and mitosis, for spindle organization and centrosome integrity. Diseases associated with CHD3 include Snijders Blok-Campeau Syndrome and Agenesis Of Corpus Callosum, Cardiac, Ocular, And Genital Syndrome.

Protein Details: Recombinant CHD3 protein was expressed in baculovirus expression system as the full length protein (accession number NP_005843.2) with a N-terminal FLAG tag. The molecular weight of the protein is 224 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 CHD3 protein

7.5% SDS-PAGE with Coomassie staining

MW: 224 kDa

Purity: >80%