

## Recombinant CHD4 protein

**Catalog No:** 81385, 81685

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.7 µg/µl

**Source:** Human

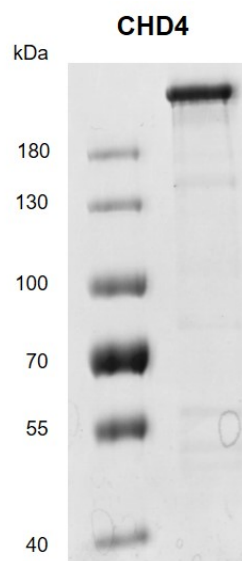
**Buffer Contents:** Recombinant CHD4 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

**Background:** CHD4 (Chromodomain Helicase DNA Binding Protein 4) 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.

**Protein Details:** Recombinant CHD4 protein was expressed in a baculovirus expression system as the full length protein (accession number NP\_001264.2) with a N-terminal FLAG tag. The molecular weight of the protein is 219.3 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 CHD4 protein**

7.5% SDS-PAGE with Coomassie staining

MW: 219.3 kDa

Purity: >90%