

PRODUCT NAME		
CHD4 polyclonal antibody		
Other names: Mi-2, Mi-2b, Mi2-BETA, CHD-4		
Cat. No. C15410031 (pAb-031-050)	Type: Polyclonal	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

Description: Polyclonal antibody raised in rabbit against human CHD4 (chromodomain helicase DNA binding protein 4), using a recombinant protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:500	Fig1

Purity: Protein G purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Last data sheet update: April 23, 2010

Target description

CHD4 (UniProtKB/Swiss-Prot entry Q14839) belongs to the SNF2/RAD54 helicase family. It is a putative transcription regulator and is a central component of the nucleosome remodeling and histone deacetylase (NuRD) repressor complex which plays an important role in epigenetic transcriptional repression. CHD4 was initially discovered as autoantigen in the connective tissue disease dermatomyositis.

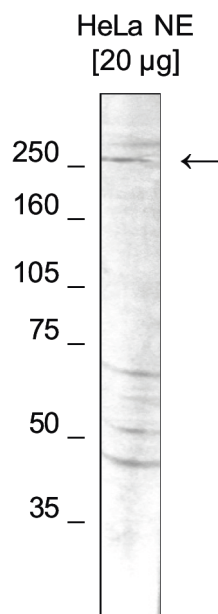


Figure 1

Western blot analysis using the Diagenode antibody directed against CHD4

Western blot was performed on nuclear extracts from HeLa cells (HeLa NE, 20 µg) using the Diagenode antibody against CHD4 (cat# pAb-031-050) diluted 1:500 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.