

PRODUCT NAME hSNF2H monoclonal antibody		
Other names: SMARCA5, ISWI, WCRF135]		
Catalog #: MAb-040-050	Type: Monoclonal IgG2b	Size: 50 µg/ 25 µl
Lot #: 001	Source: Mouse	Concentration: 2.0 µg/µl

Description: Monoclonal antibody raised in mouse against human SNF2H (SWI/SNF-related, matrix-associated, actin-dependent) using a recombinant protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1
Immunofluorescence	1:100 - 1:500	Ref 1
Immunoprecipitation	0.3 µg	Ref 1

Purity: Protein G purified monoclonal 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: February 18, 2010

References

[1] Loyola A., Huang J.Y., LeRoy G., Hu S., Wang Y.H., Donnelly R.J., Lane W.S., Lee S.C. and Reinberg D. 2003. *Mol. Cell. Biol.* 23[19]:6759-68.

Product description

Complexes containing Snf2h (SMARCA5) (UniProtKB/Swiss-Prot entry O60264) are capable of forming ordered nucleosome arrays on chromatin in vitro (may require intact histone H4 tails). SNF2h is also required for replication of pericentric heterochromatin in S-phase, specifically in conjunction with BAZ1A. Snf2h probably also plays a role in repression of polI dependent transcription of the rDNA locus, through the recruitment of the SIN3/HDAC1 corepressor complex to the rDNA promoter. Snf2h is a catalytic subunit of the four known chromatin-remodeling complexes: CHRAC, RSF, ACF/WCRF, and WICH. Each complex contains subunits which may regulate the specificity or catalytic activity of SMARCA5. Snf2h possesses intrinsic ATP-dependent nucleosome remodelling activity.

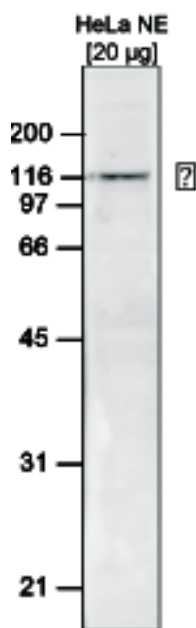


Figure 1

Western blot analysis using the Diagenode monoclonal antibody against hSnf2h

Western blot was performed on nuclear extracts from HeLa cells (HeLa NE, 20 µg) using the Diagenode monoclonal antibody directed against hSNF2H (cat# MAb-040-050), diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left, the position of the protein of interest is shown on the right.