

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

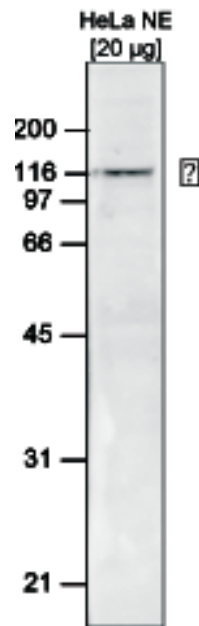
**Last data sheet update:** February 18, 2010

**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 *poll* 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.