

PRODUCT NAME		
SKI8 polyclonal antibody		
Other names: WDR61, REC14		
Cat. No. C15410012 (pAb-012-150)	Type: Polyclonal ChIP	Size: 150 µg/ 60 µl
Lot #: 001	Source: Rabbit	Concentration: 2.5 µg/µl

Description: Polyclonal antibody raised in rabbit against human SKI8, using a recombinant protein.

Specificity: Human and Drosophila: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1, Ref 1
ChIP	1 - 5 µg per ChIP	Ref 1

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.

Availability date: February 19, 2007. Last data sheet update: November 6, 2007

This antibody has been described in:

(1) Zhu B, Mandal SS, Pham AD, Zheng Y, Erdjument-Bromage H, Batra SK, Tempst P and Reinberg D (2005)
The human PAF complex coordinates transcription with events downstream of RNA synthesis. Genes Dev 19:1668-1673.

Last data sheet update: April 8, 2010

Target description

SKI8 (UniProtKB/Swiss-Prot entry Q9GZS3) is a component of the PAF and SKI complex which is composed of SKI2, SKI3 and SKI8. These are involved in transcriptional regulation and events downstream of RNA synthesis. Interaction of the SKI complex with SKI7 links the SKI complex to the exosome in order to perform mRNA degradation.

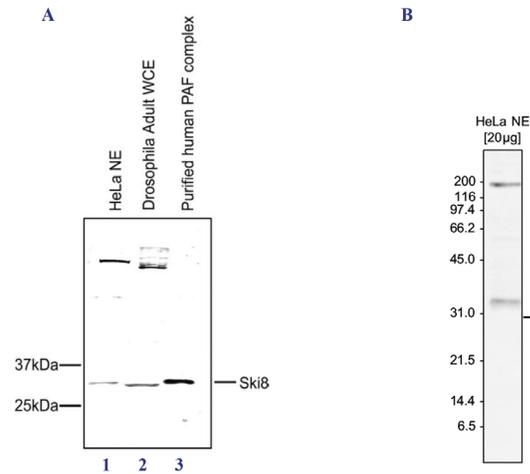


Figure 1

Western blot analysis using the Diagenode antibody directed against SKI8

Western blot was performed on nuclear extract from HeLa cells (lane 1), Drosophila adult whole cell extracts (lane 2), and the purified human PAF complex (lane 3) (Figure 1A) or on nuclear extracts from HeLa cells (HeLa NE, 20 µg) (Figure 1B), using the Diagenode antibody against SKI8 (Cat. No. pAb-012-150) diluted 1:1,000. The location of the protein is indicated on the right, the marker (in kDa) is shown on the left.