

PRODUCT NAME		
dCdc73 polyclonal antibody		
Other names: HRPT2, HPTJT, Parafibromin		
<b>Cat. No.</b> C15410019 (pAb-019-050)	<b>Type:</b> Polyclonal	<b>Size:</b> 50 µg/ 25 µl
<b>Lot #:</b> 001	<b>Source:</b> Rabbit	<b>Concentration:</b> 2.0 µg/µl

**Description:** Polyclonal antibody raised in rabbit against drosophila Cdc73 (Cell division cycle protein 73), using the full length recombinant protein.

**Specificity:** Drosophila and human: positive  
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig1
Immunofluorescence	12 µg per IP	
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.

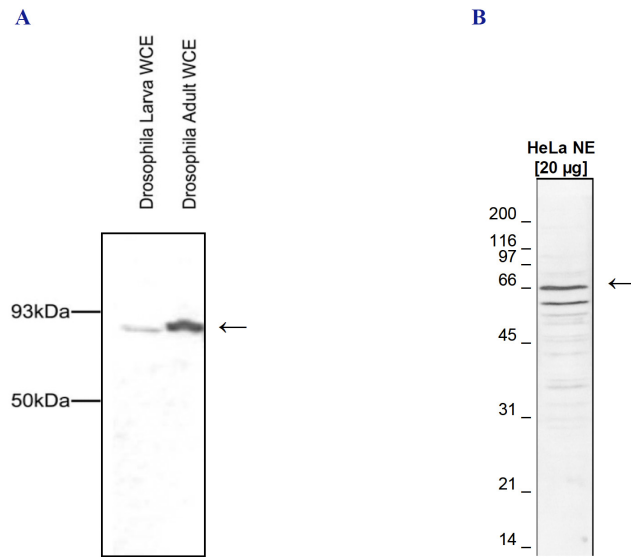
**This antibody has been described in:**

(1) Adelman K, Wei W, Ardehali MB, Werner J, Zhu B, Reinberg D and Lis JT (2006) Drosophila Paf1 modulates chromatin structure at actively transcribed genes. Mol Cell Biol 26:250-260.

**Last data sheet update:** April 9, 2010

**Target description**

Cdc73 (UniProtKB/Swiss-Prot entry Q6P1J9) is a tumor suppressor probably involved in transcriptional and post-transcriptional control pathways. Cdc73 is part of the PAF1 complex, a five-subunit protein complex composed of Paf1, Cdc73, Leo1, Rtf1 and Ctr9. PAF1 was first purified from yeast in association with RNA polymerase II and is believed to function in transcription elongation. The yeast PAF (yPAF) complex interacts with RNA polymerase II and coordinates the setting of histone marks associated with active transcription. In addition to coordinating events during transcription (initiation, promoter clearance, and elongation), hPAF also coordinates events in RNA quality control. Cdc73 may also be involved in cell cycle progression through the regulation of cyclin D1/PRAD1 expression.



**Figure 1**

**Western blot analysis using the Diagenode antibody directed against dCdc73**

Western blot was performed on whole cell extracts (WCE) from drosophila larva and drosophila adults (Figure 1A) or on nuclear extracts from HeLa cells (HeLa NE, 20 µg) (Figure 1B) using the Diagenode antibody against dCdc73 (Cat. No. pAb-019-050) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. A molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.