

| PRODUCT NAME | | |
|---|-------------------------|---------------------------------|
| dRtf1 polyclonal antibody | | |
| Other names: RDL | | |
| Cat. No. C15410018 (pAb-018-050) | Type: Polyclonal | Size: 50 µg/ 28 µl |
| Lot #: 001 | Source: Rabbit | Concentration: 1.8 µg/µl |

Description: Polyclonal antibody raised in rabbit against drosophila Rtf1 (Rtf1, Paf1/RNA polymerase II complex component, homolog), using the full length recombinant protein.

Specificity: Drosophila: positive
Other species: not tested

| Applications | Suggested dilution | References |
|---------------------|--------------------|--------------|
| Western blotting | 1:1000 | Fig 1; Ref 1 |
| Immunofluorescence | 1:50 | Ref 1 |
| Immunoprecipitation | 12 µg per IP | 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.

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

PAF is a five-subunit protein complex composed of Paf1, Cdc73, Leo1, Rtf1 and Ctr9, which was first purified from yeast. The yeast PAF (yPAF) complex interacts with RNA polymerase II and coordinates the setting of histone marks associated with active transcription. The human PAF (hPAF) complex shares four subunits with yPAF (hCtr9, hPaf1, hLeo1 and Cdc73), but contains an additional higher eukaryotic-specific subunit, hSki8. In addition to coordinating events during transcription (initiation, promoter clearance and elongation); PAF also coordinates events in RNA quality control.



Figure 1

Western blot analysis using the Diagenode antibody directed against dRtf1

Whole cell extracts (WCE) from drosophila larva and drosophila adults were analysed by Western blot using the Diagenode antibody against dRtf1 (Cat. No. pAb-018-050), diluted 1:1000. Size and location of the protein are indicated.