

PRODUCT NAME SPT16 polyclonal antibody		
Other names: SUPT16H, FACT140, CDC68, FACTP140		
Cat. No. C15410049 (pAb-049-050)	Type: Polyclonal	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

Description: Polyclonal antibody raised in rabbit against human SPT16 (Suppressor of Ty 16 homolog), using a recombinant protein containing amino-acids 276 to 572.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1; [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] Orphanides G, Wu WH, Lane WS, Hampsey M and Reinberg D (1999) The chromatin-specific transcription elongation factor FACT comprises human SPT16 and SSRP1 proteins. Nature 400: 284-288.

Last data sheet update: April 21, 2010

Target description

SPT16 (UniProtKB/Swiss-Prot entry Q9Y5B9) is a component of the FACT complex, a general chromatin factor that acts to reorganize nucleosomes. The FACT complex is involved in multiple processes that require DNA as a template such as mRNA elongation, DNA replication and DNA repair. During transcription elongation the FACT complex acts as a histone chaperone that both destabilizes and restores nucleosomal structure. SPT16 facilitates the passage of RNA polymerase II and transcription by promoting the dissociation of one histone H2A-H2B dimer from the nucleosome and subsequently promotes the reestablishment of the nucleosome. The FACT complex is probably also involved in phosphorylation of 'Ser-392' of p53/TP53 via its association with casein kinase II.

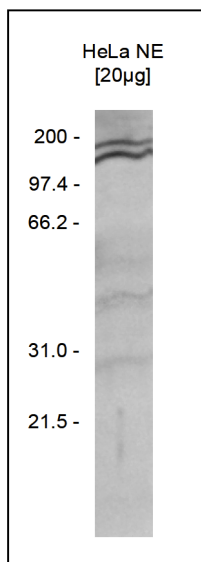


Figure 1

Western blot analysis using the Diagenode antibody directed against SPT16

HeLa nuclear extracts (HeLa NE, 20 µg) were analysed by Western blot using the Diagenode antibody against SPT16 (Cat. No. pAb-049-050) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. A molecular weight marker is shown on the left.