

PRODUCT NAME		
SSRP1 polyclonal antibody		
Other names: FACT80, FACTp80		
Cat. No. C15410050 (pAb-050-050)	Type: Polyclonal	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

Description: Polyclonal antibody raised in rabbit against human SSRP1 (Structure-specific recognition protein 1), the GST tagged N-terminal part of the protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 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.

Last data sheet update: March 17, 2010

Target description

SSRP1 (UniProtKB/Swiss-Prot entry Q08945) is a component of the FACT complex, a general chromatin factor that acts to reorganize nucleosomes. The FACT complex, which besides SSRP1 also contains SUPT16H (SPT16) 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. It facilitates the passage of RNA polymerase II and transcription by promoting the dissociation of one histone H2A-H2B dimer from the nucleosome, subsequently promoting the reestablishment of the nucleosome following the passage of RNA polymerase II. The FACT complex is probably also involved in phosphorylation of 'Ser-392' of p53/TP53 via its association with CK2 (casein kinase II). SSRP1 binds specifically to double-stranded DNA and at low levels to DNA modified by the antitumor agent cisplatin. SSRP1 may potentiate cisplatin-induced cell death by blocking replication and repair of modified DNA. SSRP1 also acts as a transcriptional coactivator for p63/TP73L.

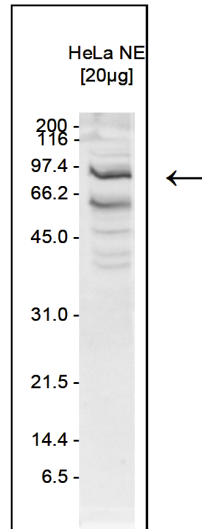


Figure 1

Western blot analysis using the Diagenode antibody directed against hSSRP1

Western blot was performed on nuclear extracts from HeLa cells (20 µg) using the Diagenode antibody against hSSRP1 (Cat. No. pAb-050-050) 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 by an arrow.