

PRODUCT NAME		
Set9 polyclonal antibody		
Other names: SET7, SET7/9, KMT7, H3-K4-HMTase		
Catalog #: C15310047 (CS-047-100)	Type: Polyclonal	Size: 100 µl
Lot #: 001	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human Set9 (SET domain containing (lysine methyltransferase) 9), using a recombinant protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig1; [1]

Purity: Whole antiserum from rabbit containing 0.05% azide.

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: April 12, 2010

Target description

Set9 (UniProtKB/Swiss-Prot entry Q8WTS6) is a histone methyltransferase which specifically methylates lysine 4 of histone H3 being a specific tag for epigenetic transcriptional activation. Further, methylation of histone H3K4 by Set9 precludes deacetylation of H3 by the NuRD complex and impairs Suv39h1-mediated methylation of H3K9. Set9 also methylates the tumor suppressor p53 at lysine 372, thereby enhancing its stability. It contains a SET domain but lacks the pre- and post-SET domains.).

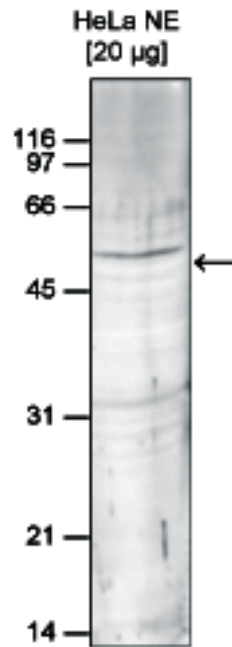


Figure 1

Western blot analysis using the Diagenode antibody directed against Set9

Western blot was performed on nuclear extracts from HeLa cells (HeLa NE, 20 µg) using the Diagenode antibody against Set9 (cat# CS-047-100) 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.