

## SETDB1 polyclonal antibody - Classic

**Other names:** ESET, KMT1E, KG1T, TDRD21, H3-K9-HMTase4

**Cat. No.** C15410249

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 40471

**Size:** 25 µl/100 µl

**Concentration:** 0.63 µg/µl

**Specificity:** Human.

**Purity:** Purified by antigen-affinity chromatography.

**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.

### Applications

	Suggested dilution*	Results
Western blot	1:500-1:3,000	Figure 1
Immunoprecipitation	1:100-1:500	Figure 2

\*Optimal dilutions/concentrations should be determined by the researcher.

### Target description

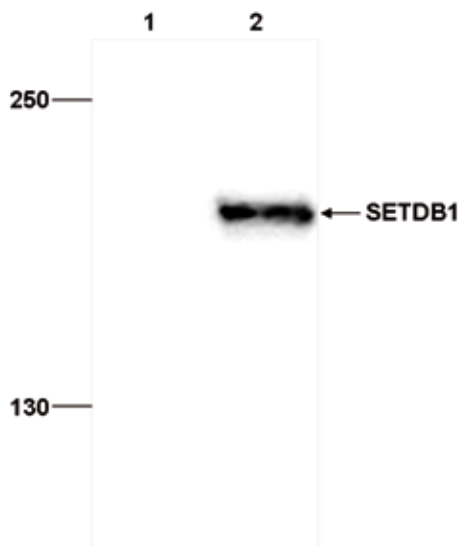
SETDB1 (UniProt/Swiss-Prot entry Q15047) is a Histone methyltransferase that specifically trimethylates Lys9 of histone H3 thereby playing a central role in transcriptional repression and euchromatin formation. H3K9 trimethylation is coordinated with DNA methylation, probably by association of SETDB1 with MBD1 and ATF7IP. SETDB1 has been identified as a target for treatment in Huntington Disease since gene silencing and transcription dysfunction likely play a role in the disease pathogenesis.

Results



**Figure 1. Western Blot analysis using the Diagenode antibody directed against SETDB1**

Whole cell extracts (30 µg) from HeLa cells were analysed by Western blot using the Diagenode antibody against SETDB1 (Cat. No. C15410249) diluted 1:10,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 2. Immunoprecipitation using the Diagenode antibody directed against SETDB1**

Immunoprecipitation was performed on whole cell extracts from HeLa cells using 4 µg of the Diagenode antibody against SETDB1 (Cat. No. C15410249, lane 1). An equal amount of rabbit IgG was used as a negative control (lane 2). The immunoprecipitated SETDB1 protein was detected by western blot with the SETDB1 antibody diluted 1:1,000.