

PRODUCT NAME		
Setd1b polyclonal antibody		
Other names: SET1B, KMT2G		
<b>Cat. No.</b> C15310118 (CS-118-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A278-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse Setd1b (Set domain containing 1B), using a 3 KLH-conjugated synthetic peptides containing sequences from the central part of the protein (1).

**Specificity:** Mouse: positive  
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:500	Fig 2 (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.

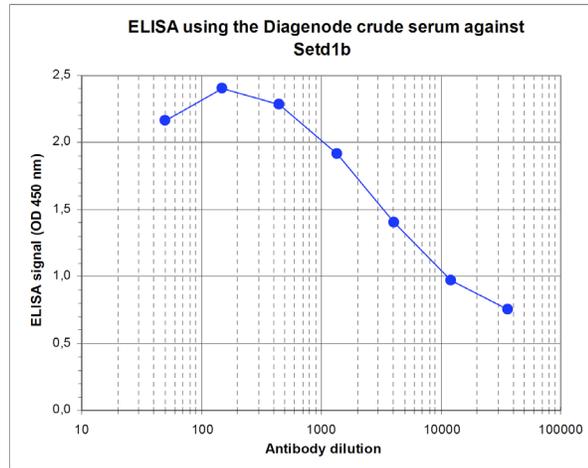
**References:**

(1) Peptide design by Andrea Kranz, Western blot analysis by Heike Petzold and Andrea Kranz, BIOTEC, Dept. of Genomics, Prof. A. F. Stewart, TU Dresden, Tatzberg 47/49, 01307 Dresden, Germany

**Last data sheet update:** March 23, 2010

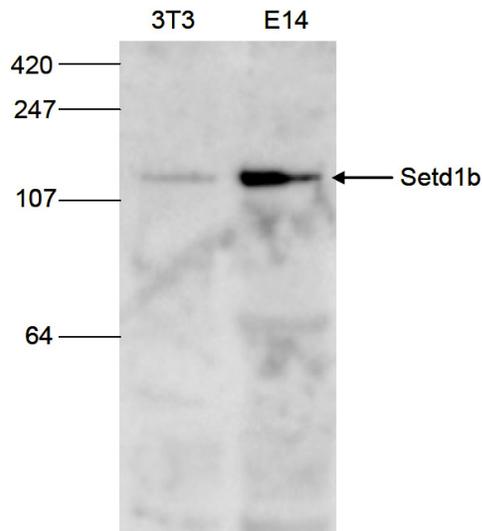
**Target description**

SETD1B (UniProt/Swiss-Prot entry Q9UPS6) is a component of the SET1 histone methyltransferase (HMT) complex. This complex specifically methylates lysine 4 of histone H3, but only if the neighboring lysine 9 residue is not yet methylated. Methylation of H3K4 represents a specific tag for epigenetic transcriptional activation. SETD1B shows a non-overlapping localization with SETD1A, suggesting that both proteins play a specific role in the epigenetic control of chromatin structure and gene expression.



**Figure 1**  
**Determination of the titer**

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse Setd1b [Cat. No. CS-118-100]. The plates were coated with a mix of the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:8,400.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against mouse Setd1b (1)**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Setd1b [Cat. No. CS-118-100], diluted 1:500 in BSA/PBS-Tween. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.