

PRODUCT NAME		
<b>Dnmt2 polyclonal antibody</b>		
Other names: TRDMT1, M.HsaIIP, PuMet, RNMT1		
<b>Cat. No.</b> C15310095 (CS-095-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A359-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse Dnmt2 (DNA methyltransferase 2), using a KLH-conjugated synthetic peptide containing an amino acid sequence from the N-terminal part of the protein [1].

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

Applications	Suggested dilution	References
ELISA	1:100 – 1:1,000	Fig 1
Western blotting	1:1,000	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.

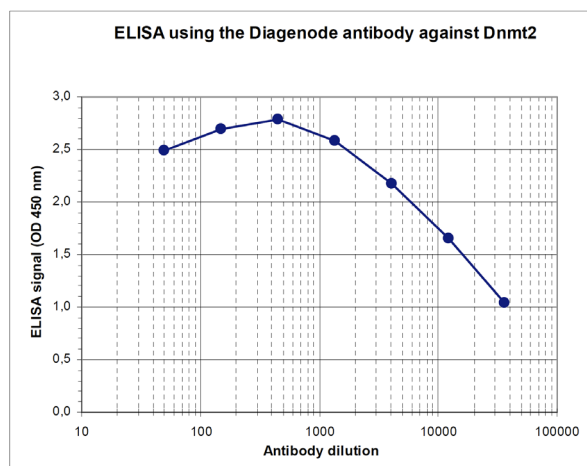
**References:**

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

**Last data sheet update:** April 7, 2010

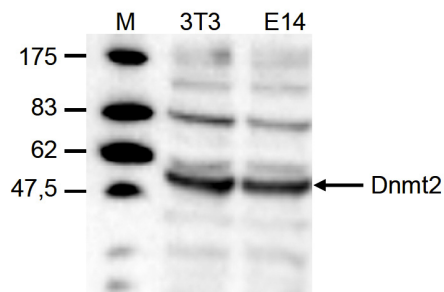
**Target description**

Dnmt2 [UniProtKB/Swiss-Prot entry O14717] belongs to the family of DNA cytosine methyltransferases. Dnmt2, however, lacks the large N-terminal regulatory domain common to other eukaryotic methyltransferases and does not methylate CpG in genomic DNA. Instead, it specifically methylates cytosine 38 in the anticodon loop of Asp transfer RNA.



**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 Dnmt2 (Cat. No. CS-095-100). The wells were coated with the peptide used for immunisation of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:21,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against Dnmt2 [1]**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Dnmt2 (Cat. No. CS-095-100), diluted 1:1000 in BSA/PBS-Tween. The molecular weight marker (M, in kDa) is shown on the left; the location of the protein of interest (47 kDa) is indicated on the right.