

PRODUCT NAME		
MLL2 polyclonal antibody		
Other names: TRX2, WBP7, MLL4		
<b>Cat. No.</b> C15310099 (CS-099-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A268-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse MLL2 (Myeloid/lymphoid or mixed-lineage leukemia protein 2), using two KLH-conjugated synthetic peptides containing an amino acid sequence from the central and from the N-terminal part of the protein, respectively (1).

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

Applications	Suggested dilution	References
ELISA	1:200 – 1:2,000	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.

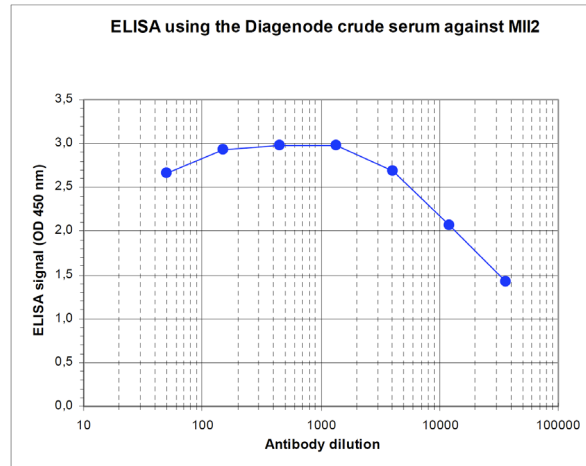
**References:**

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

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

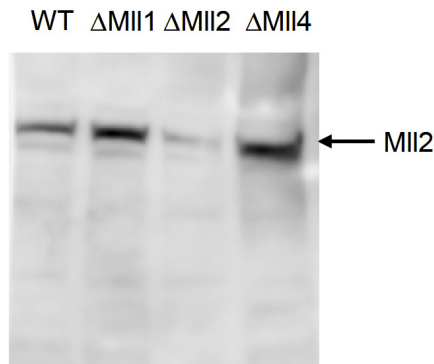
**Target description**

MLL2 (UniProtKB/Swiss-Prot entry Q9UMN6) plays a role in H3K4 methylation and may therefore act as a transcriptional regulator. MLL2 is essential for embryonic development and may also be important for hematopoiesis. Further, MLL2 shows an elevated expression in solid tumor cell lines, and may be involved in cancer.



**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 MII2 (Cat. No. CS-099-100). The wells were coated with the peptides 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 1:32,500.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against MII2 (1)**

Western blot was performed on whole cell lysates from mouse embryonic stem cells [E14Tg2a] with the Diagenode antibody against mouse MII2 (Cat. No. CS-099-100), diluted 1:500 in BSA/PBS-Tween. The location of the N-terminal fragment of MII2 is indicated on the right. Cells homozygous for the targeted conditional mII2 allele ( $\Delta$ MII2) show a dramatic reduction of MII2 protein after partial recombination, whereas MII2 protein was detected in  $\Delta$ MII1,  $\Delta$ MII4 and WT E14Tg2a cells.