

PRODUCT NAME		
MLL4 polyclonal antibody		
Other names: ALR, MLL2		
<b>Cat. No.</b> C15310100 (CS-100-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A274-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse MLL4 (Myeloid/lymphoid or mixed-lineage leukemia protein 4), using three KLH-conjugated synthetic peptides containing an amino acid sequence 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.

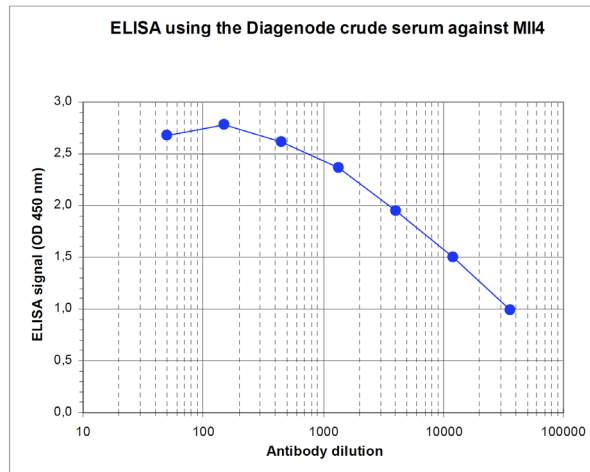
**References:**

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

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

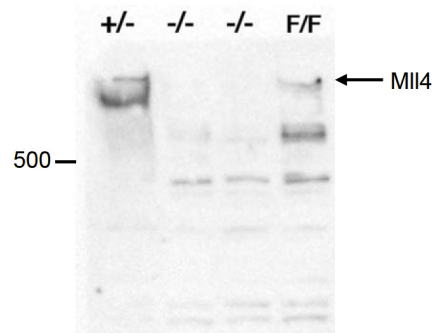
**Target description**

MLL4 (UniProtKB/Swiss-Prot entry O14686) is involved in the methylation of lys 4 of histone H3, thus acting as a transcriptional regulator.



**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 Mll4 (Cat. No. CS-100-100). 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 1:15,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against Mll4 (1)**

Western blot was performed on whole cell lysates from mouse embryonic stem cells [E14Tg2a] with the Diagenode antibody against mouse Mll4 (Cat. No. CS-100-100), diluted 1:500 in PBS-Tween containing 5% skimmed milk. The location of the protein of interest (~ 600 kDa) is indicated on the right; the location of the 500 kDa MW marker is shown on the left. Cells homozygous for the targeted mll4 allele (-/-) show a complete loss of Mll4 protein, whereas the Mll4 protein was detected in flp-recombined ES cells (F/F) and heterozygously targeted ES cells (+/-).