

MLL2 antibody (other name: TRX2, WBP7, MLL4)

Cat. No. C15310099

Type: Polyclonal

Source: Rabbit

Lot: A268-004

Size: 100 µl

Concentration: Not determined

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

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.

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 N-terminal and one peptide containing a sequence from the central part of the protein, respectively.

Applications

Applications	Suggested dilution	References
ELISA	1:200 - 1:2,000	Fig 1
Western blotting **	1:500	Fig 2, [1]

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.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: August, 2022

Results

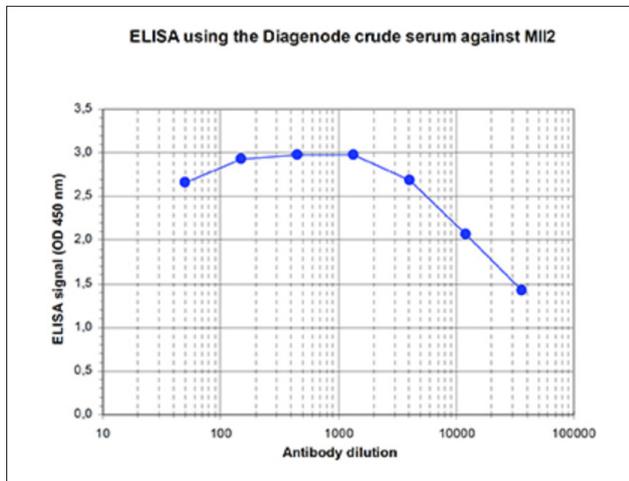


Figure 1. Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse Mll2 (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.

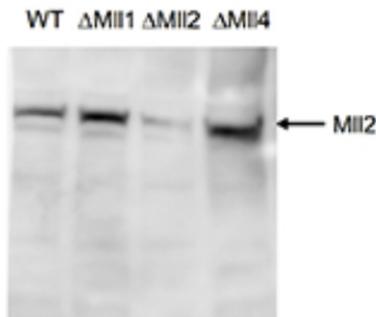


Figure 2. Western blot analysis using the Diagenode antibody directed against Mll2

Western blot was performed on whole cell lysates from mouse embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Mll2 (Cat. No. CS-099-100), diluted 1:500 in BSA/PBS-Tween. The location of the N-terminal fragment of Mll2 is indicated on the right (270 kDa). Cells homozygous for the targeted conditional mll2 allele (ΔMll2) show a dramatic reduction of Mll2 protein after partial recombination, whereas Mll2 protein was detected in ΔMll1, ΔMll4 and WT E14Tg2a cells.