

MII2 polyclonal antibody - Pioneer

Cat. No. C15310099

Type: Polyclonal	Specificity: Mouse: positive. Other species: not tested.
Size: 100 µl	Isotype: NA
Concentration: Not determined	Source: Rabbit
Lot No.: A268-004	Purity: Whole antiserum from rabbit containing 0.05% azide.
Storage buffer: NA	Storage conditions: 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

Alternative names: **TRX2, WBP7, MLL4**

Polyclonal antibody raised in rabbit against mouse MII2 (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.

Applications

Applications	Suggested dilution	References
ELISA	1:200 - 1:2,000	Fig 1
Western Blotting	1:500	Fig 2, (1)

Target Description

MII2 (UniProtKB/Swiss-Prot entry Q9UMN6) plays a role in H3K4 methylation and may therefore act as a transcriptional regulator. MII2 is essential for embryonic development and may also be important for hematopoiesis. Further, MII2 shows an elevated expression in solid tumor cell lines, and may be involved in cancer.

Validation Data

ELISA using the Diagenode crude serum against MII2

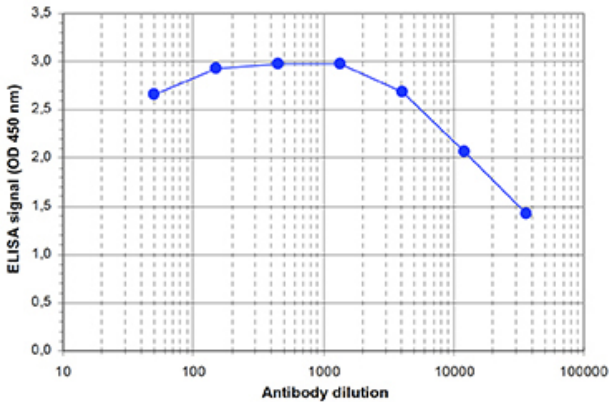


Figure 1. Determination of the titer

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.

WT ΔMII1 ΔMII2 ΔMII4

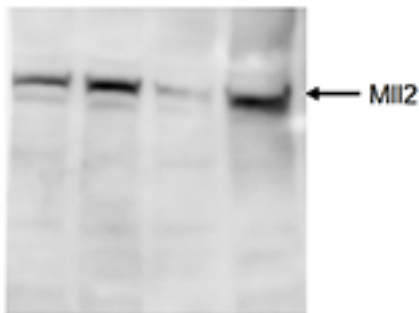


Figure 2. Western blot analysis using the Diagenode antibody directed against MII2

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 (270 kDa). Cells homozygous for the targeted conditional mii2 allele (?MII2) show a dramatic reduction of MII2 protein after partial recombination, whereas MII2 protein was detected in ?MII1, ?MII4 and WT E14Tg2a cells.