

PRODUCT NAME		
Brd2 polyclonal antibody		
Other names: FSRG1, NAT, RING3, RNF3		
Cat. No. CS-094-100	Type: Polyclonal	Size: 100 µl
Lot #: A355-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against mouse Brd2 (bromodomain containing 2), using a KLH-conjugated synthetic peptide 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	100	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

Brd2 [UniProtKB/Swiss-Prot entry P25440], a mitogen-activated kinase localized to the nucleus, is a putative transcriptional regulator which interacts with E2F1 and with histone H4 acetylated at 'Lys-13'. It is expressed during development and may be involved in growth control. Brd2 also may play a role in spermatogenesis or folliculogenesis and may be involved in some types of leukemia and in juvenile myoclonic epilepsy.

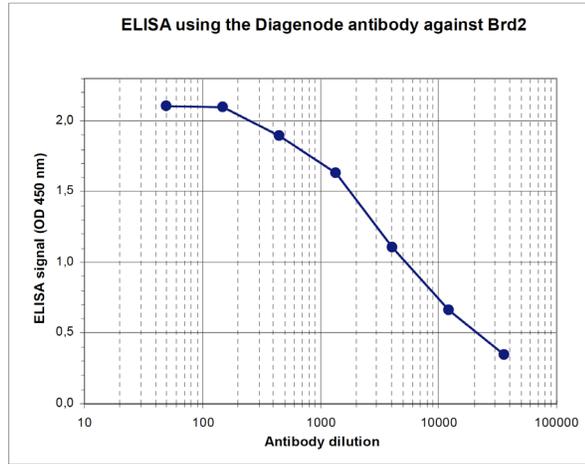


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse Brd2 (Cat. No. CS-094-100). 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 1:5,000.

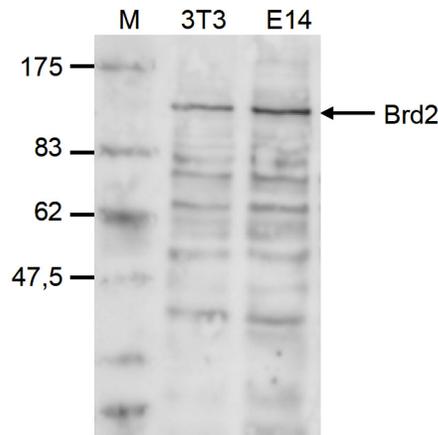


Figure 2
Western blot analysis using the Diagenode antibody directed against Brd2

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Brd2 (Cat. No. CS-094-100), diluted 1:1,000 in BSA/PBS-Tween. The molecular weight marker (M, in kDa) is shown on the left; the location of the protein of interest (predicted size: 88 kDa) is indicated on the right.