

PRODUCT NAME		
ZMYND8 polyclonal antibody		
Other names: PRKCBP1, RACK7, PR02893		
Cat. No. C15310136 (CS-136-100)	Type: Polyclonal ChIP-grade	Size: 100 µl
Lot #: A331-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human ZMYND8 (zinc finger, MYND-type containing 8), using a KLH-conjugated synthetic peptide containing a sequence from the C-terminal part of the protein.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:1,000	Fig 2

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.

Last data sheet update: March 18, 2010

Target description

ZMYND8 (UniProtKB/Swiss-Prot entry Q9ULU4) is a receptor for activated protein kinase C (RACK). It also contains a bromodomain and two zinc fingers, and is thought to be a transcriptional regulator. Further, ZMYND8 is a cutaneous T-cell lymphoma-associated antigen.

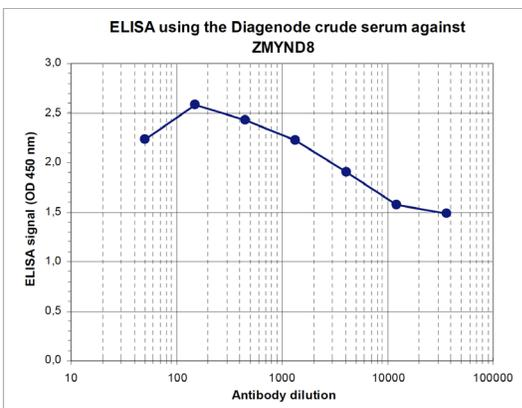


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human ZMYND8 (Cat. No. CS-136-100). The plates were coated with the peptide used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:36,000.

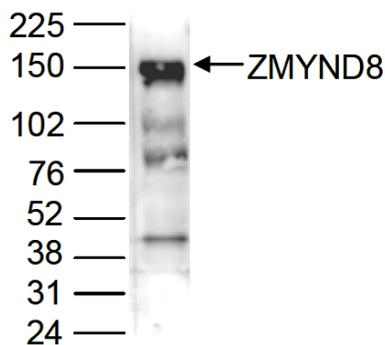


Figure 2
Western blot analysis using the Diagenode antibody directed against ZMYND8

Nuclear extracts from HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against ZMYND8 (Cat. No. CS-136-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right (expected size: 132 kDa); the marker (in kDa) is shown on the left.