

PRODUCT NAME		
ZBTB38 polyclonal antibody		
Other names: CIBZ		
Cat. No. C15310123 (CS-123-100)	Type: Polyclonal	Size: 100 µl
Lot #: A233-001	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human ZBTB38 (zinc finger and BTB domain containing 38), using three KLH-conjugated synthetic peptides, two containing a sequence from the central part and one 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 – 1:1,000	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 22, 2010

Target description

ZBTB38 (UniProt/Swiss-Prot entry Q8NAP3) is a methyl CpG binding protein that binds to DNA sequences containing a single methylated CpG, thereby acting as a transcriptional repressor. It has been found to be highly expressed in the brain and may be involved in the differentiation of late postmitotic neurons.

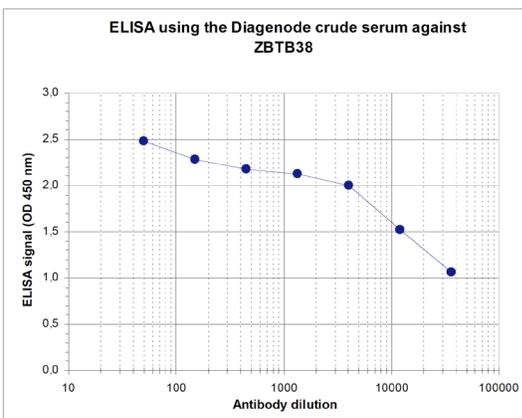


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human ZBTB38 [Cat. No. CS-123-100]. The plates were coated with the peptides 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 1:26,000.

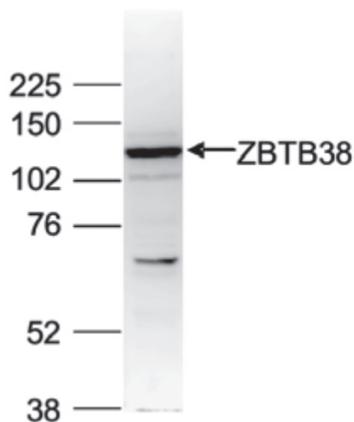


Figure 2
Western blot analysis using the Diagenode antibody directed against ZBTB38

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against ZBTB38 [Cat. No. CS-123-100] diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest (expected size: 138 kDa) is indicated on the right; the marker (in kDa) is shown on the left.