

| PRODUCT NAME                           |   |                                      |
|--|---|--------------------------------------|
| ZBTB4 polyclonal antibody              |   |                                      |
| Other names: KAISO-L1                  |   |                                      |
| <b>Cat. No.</b> C15310142 (CS-142-100) | <b>Type:</b> Polyclonal <b>ChIP-grade</b> | <b>Size:</b> 100 µl                  |
| <b>Lot #:</b> A452-001                 | <b>Source:</b> Rabbit                     | <b>Concentration:</b> not determined |

**Description:** Polyclonal antibody raised in rabbit against human ZBTB4 (zinc finger and BTB domain containing 4), using four different KLH-conjugated synthetic peptides.

**Specificity:** Human: positive  
Other species: not tested

| Applications     | Suggested dilution | References |
|------------------|--------------------|------------|
| ELISA            | 1:1,000 – 1:5,000  | Fig 1      |
| Western blotting | 1:1,000            | Fig 2      |

**Purity:** Whole antiserum from rabbit containing 0.05% azide.

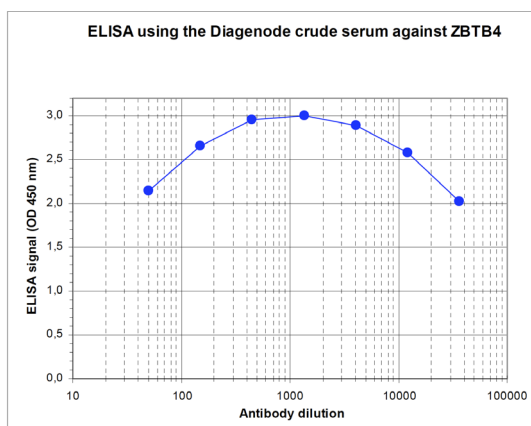
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** March 3, 2010

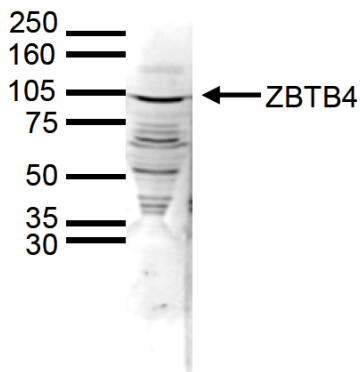
**Target description**

ZBTB4 (UniProtKB/Swiss-Prot entry Q9P1Z0) is a KAISO like protein which binds to methylated DNA and regulates the transcription of methylated targets. Unlike KAISO, ZBTB4 is able to bind to single methylated CpG's. ZBTB4 acts as a transcriptional repressor protein and has been shown to be downregulated in the advanced stage of different tumours.



**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 ZBTB4 (Cat. No. CS-142-100). 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 1:50,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against ZBTB4**

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against ZBTB4 (Cat. No. CS-142-100) 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: 105 kDa); the marker (in kDa) is shown on the left.