

## CBX3 polyclonal antibody

**Other names:** HP1gamma, HECH

**Cat. No.** C15310234 (CS-PA014-100)

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** A1131-001

**Size:** 100 µl

**Concentration:** not determined

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

**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.

**Description:** Polyclonal antibody raised in rabbit against human CBX3 (chromobox homolog 3) using two KLH-conjugated synthetic peptides containing a sequence from the N-terminal and the central region of the protein, respectively.

### Applications

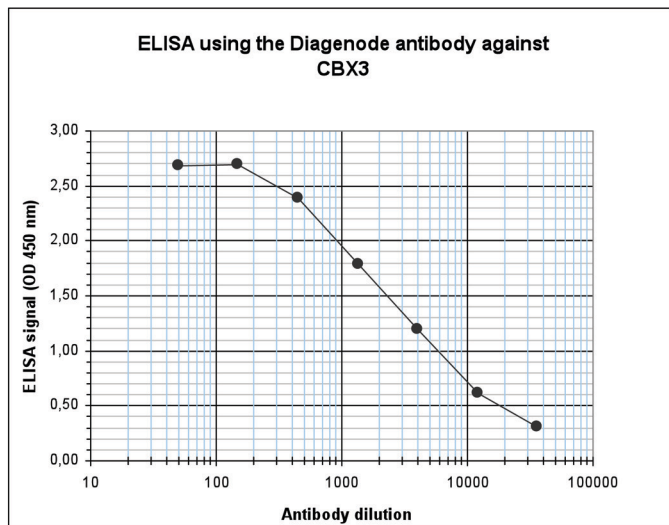
	Suggested dilution	Results
ELISA	1:100	Fig 1

\*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:500 dilution

### Target description

CBX3 (UniProtKB/Swiss-Prot entry Q13185) is involved in transcriptional silencing in heterochromatin-like complexes. Epigenetic repression is established by binding of CBX3 to histone H3 methylated at lysine 9. CBX3 may contribute to the association of the heterochromatin with the inner nuclear membrane through its interaction with lamin B receptor (LBR) and seems to be involved in the formation of functional kinetochore through interaction with MIS12 complex.

## Results



**Figure 1. Determination of the antibody titer**

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human CBX3 (cat. No. CS-PA014-100). 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 1:3,250.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** September 4, 2012