

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

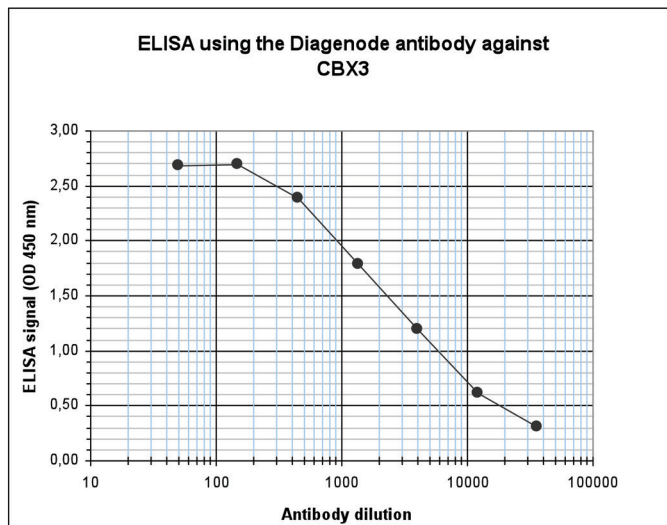


Figure 1. Determination of the antibody titer

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