

PRODUCT NAME		
HP1 alpha, beta and gamma polyclonal antibody		
Other names: CBX5, 1 and 3		
Cat. No. C15310071 (CS-071-100)	Type: Polyclonal	Size: 100 µl
Lot #: 001	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human HP1 beta (Heterochromatin protein 1 homolog beta), using the full length recombinant GST tagged protein. The antibody also recognizes the alpha and gamma isoforms.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1

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: April 13, 2010

Target description

HP1 alpha, beta and gamma (UniProtKB/Swiss-Prot entry P45973, P83916 and Q13185) are components of heterochromatin. They recognize and bind histone H3 tails methylated at 'Lys-9', leading to epigenetic repression. HP1 alpha, beta and gamma also interact with lamin B receptor (LBR) and can contribute to the association of the heterochromatin with the inner nuclear membrane.

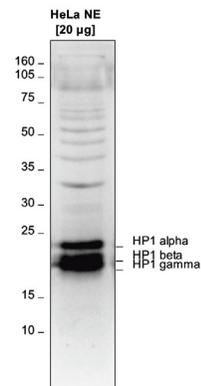


Figure 1

Western blot analysis using the Diagenode antibody against hHP1 alpha, beta and gamma

Western blot was performed on nuclear extracts from HeLa cells (HeLa NE, 20 µg) using the Diagenode antibody against HP1 alpha, beta and gamma (Cat. No. CS-071-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. A molecular weight marker (in kDa) is shown on the left. The location of the three HP1 isoforms is indicated on the right.