

PRODUCT NAME		
Suz12 pyclonal antibody		
Other names: CHET9, JJAZ1		
Cat. No. C15410029 (pab-029-050)	Type: Polyclonal ChIP-grade	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

Product description: Polyclonal antibody raised in rabbit against the N-terminus of recombinant human Suz12 protein (Suppressor of Zeste 12 protein homolog).

Specificity: Human and mouse: positive
Other species: not tested

Applications	Suggested dilution	References
ChIP*	10 µg/ChIP	Fig1
Western blotting	1:1,000	Fig 2

*Please note that of the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-10 µg per IP.

Purity: Protein G purified polyclonal antibody in PBS including 0.05% azide and 0.05% ProClin 300.

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 28, 2011

Target description

Suz12 (UniProtKB/Swiss-Prot entry Q15022) is a Polycomb group (PcG) protein. These proteins form multiprotein complexes, which keep homeotic genes in a transcriptionally repressive state throughout development. PcG proteins are not required to initiate repression, but maintain it during later stages of development. They probably act via the methylation of histones. Suz12 is a component of the PRC2 complex, which is composed of EED, EZH2, SUZ12/JJAZ1, RBBP4 and RBBP7, and methylates 'Lys-9' and 'Lys-27' of histone H3.

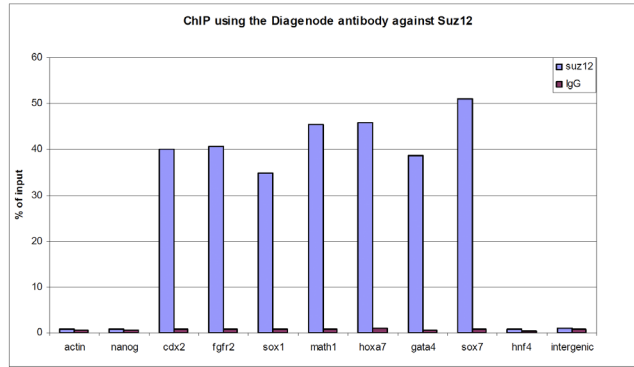


Figure 1

ChIP using the Diagenode antibody directed against Suz12

ChIP was performed using 600 µg of chromatin from mouse ES cells and 10 µg of the Diagenode antibody against Suz12 (Cat. No. pAb-029-050). IgG (10 µg) was used as a negative IP control. The IP'd DNA was analysed by QPCR with specific primer sets for different positive and negative genes. Figure 1 shows the results expressed as a % of input (the relative abundance of target DNA in the immunoprecipitates compared to the input DNA).

Data provided by A. Helness, Epigenetics and Development Group, Faculty of Medicine Imperial College London, UK.

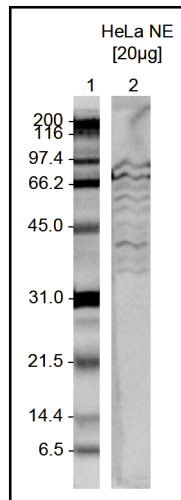


Figure 2

Western blot analysis using the Diagenode antibody directed against Suz12

HeLa nuclear extracts (HeLa NE, 20 µg) were analysed by Western blot using the Diagenode antibody directed against Suz12 (Cat. No. pAb-029-050) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk (lane 2, expected size: 83 kDa). A molecular weight marker is shown in lane 1.