

SAP30 polyclonal antibody - Classic

Cat. No. C15410036

Type: Polyclonal ChIP-grade/ChIP-seq grade

Source: Rabbit

Lot #: 001

Size: 50 µl/25 µl

Concentration: 2.0 µg/µl

Specificity: Human: positive

Other species: not tested

Purity: Protein G purified polyclonal antibody in PBS containing 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.

Description: Polyclonal antibody raised in rabbit against human SAP30 (Sin3-associated polypeptide, 30kDa), using the full length His-tagged protein.

Applications

	Suggested dilution	References
ChIP*	2 - 5 µg/ChIP	Fig 1, 2
Western blotting	1:1,000	Fig 3

*Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 µg per IP.

Target description

SAP30 (UniProtKB/Swiss-Prot entry O75446) is a component of the histone deacetylase complex, which also includes SIN3, SAP18, HDAC1, HDAC2, RbAp46, RbAp48. Histone acetylation and deacetylation play a key role in the regulation of eukaryotic gene expression. This complex only deacetylates core histone octamers, and is not able to deacetylate nucleosomal histones.

Results

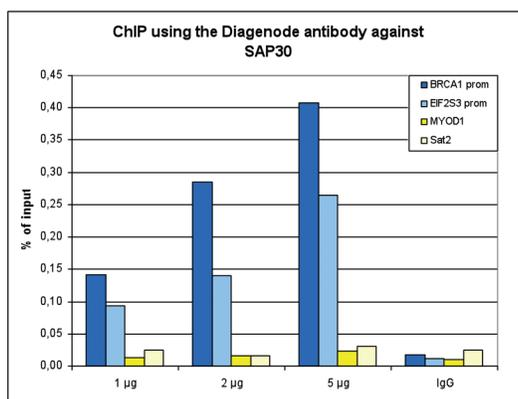


Figure 1. ChIP results obtained with the Diagenode antibody directed against SAP30

ChIP assays were performed using HeLa cells, the Diagenode antibody against SAP30 (Cat. No. C15410036) and optimized primer sets for qPCR. ChIP was performed with the "iDeal ChIP-seq" kit (cat. No. C01010055), using sheared chromatin from 4 million cells. A titration of the antibody consisting of 1, 2 and 5 µg per ChIP experiment was analysed. IgG (1 µg/IP) was used as negative IP control. QPCR was performed with primers for the EIF2S3 and BRCA1 promoters, used as positive controls, and for the MYOD1 gene and the Sat2 satellite repeat, used as negative controls. Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).

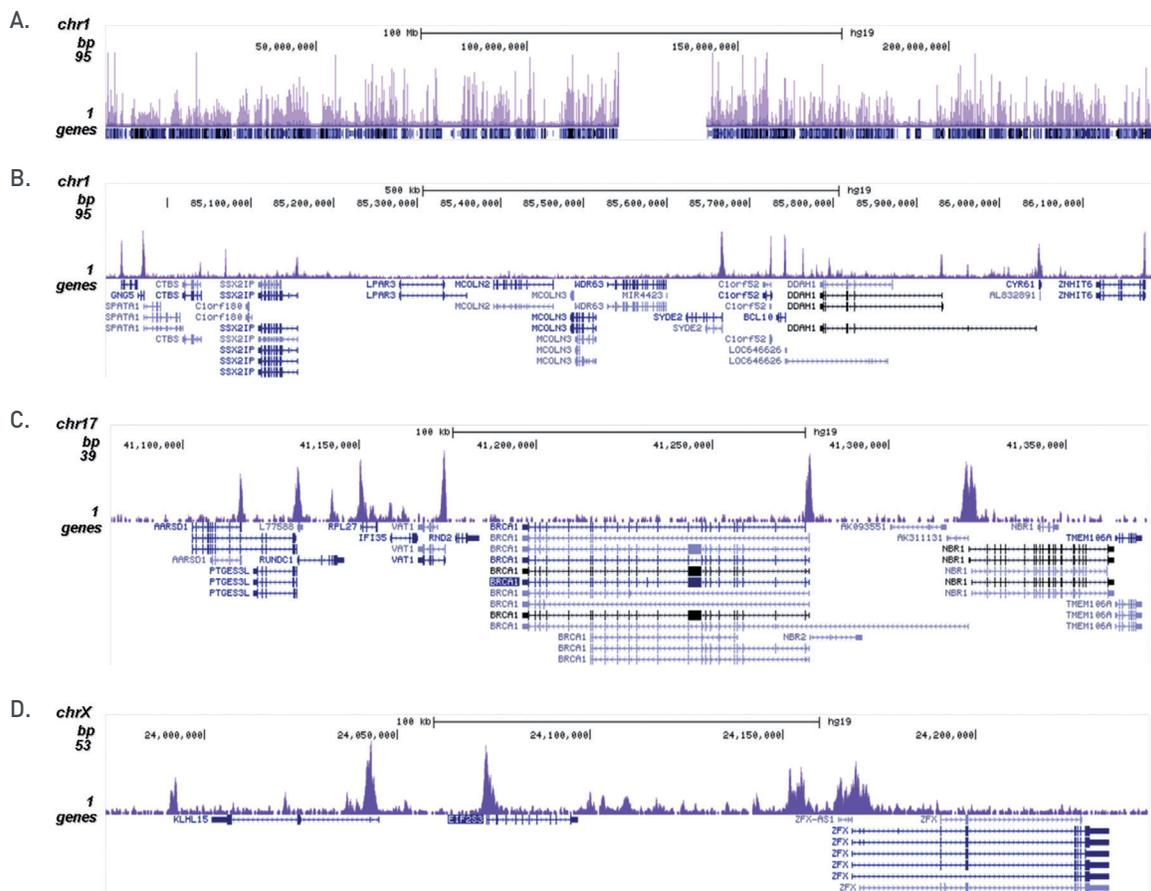


Figure 2. ChIP-seq results obtained with the Diagenode antibody directed against SAP30

ChIP was performed on sheared chromatin from 4 million HeLa cells as described above using 5 µg of the Diagenode antibody against SAP30 (Cat. No. C15410036). The IP'd DNA was subsequently analysed on an Illumina HiSeq. Library preparation, cluster generation and sequencing were performed according to the manufacturer's instructions. The 50 bp tags were aligned to the human genome using the BWA algorithm. Figure 2 shows the enrichment along the complete sequence and a 1.5 Mb region of human chromosome 1 (fig 2A and B) and in two genomic regions surrounding the BRCA1 and EIF2S3 genes on chromosome 17 and X, respectively (fig 2C and D).

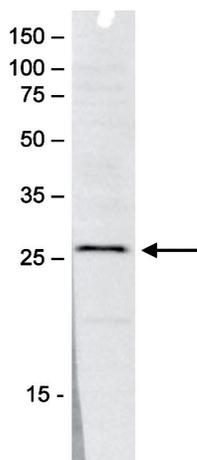


Figure 3. Western blot analysis using the Diagenode antibody directed against SAP30

Western blot was performed on nuclear extracts from HeLa cells (20 µg) using the Diagenode antibody against SAP30 (Cat. No. C15410036) diluted 1:1000 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest is indicated on the right.

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: october 3, 2015