

## p300 antibody (mAb)

**Catalog Nos:** 61401, 61903, 61402

**RRID:** AB\_2716754

**Clone:** NM11

**Isotype:** IgG1

**Application(s):** ChIP, ChIP-Seq, ICC, IF, IP, WB

**Reactivity:** Human

**Quantities:** 100 µg, 50 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 300 kDa

**Background:** p300 (E1A binding protein p300) functions as a histone acetyltransferase and regulates transcription via chromatin remodeling. Acetylates all four core histones in nucleosomes. Histone acetylation gives an epigenetic tag for transcriptional activation. Mediates cAMP-gene regulation by binding specifically to phosphorylated CREB protein. Also functions as an acetyltransferase for nonhistone targets. Acetylates 'Lys-131' of ALX1 and acts as its coactivator in the presence of CREBBP. Acetylates SIRT2 and is proposed to indirectly increase the transcriptional activity of TP53 through acetylation and subsequent attenuation of SIRT2 deacetylase function. Acetylates HDAC1 leading to its inactivation and modulation of transcription. Can also mediate transcriptional repression. Binds to and may be involved in the transforming capacity of the adenovirus E1A protein.

**Immunogen:** The antibody was raised against full-length p300 purified from human 293 cells.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

ChIP: 5 µg per ChIP

ChIP-Seq: 5 µg each

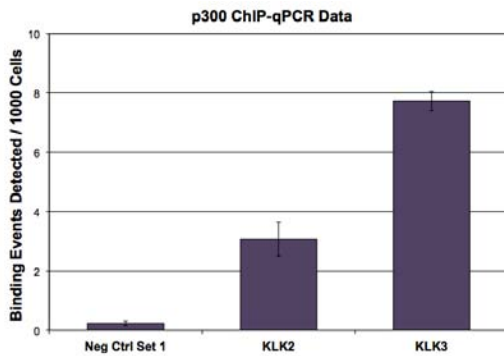
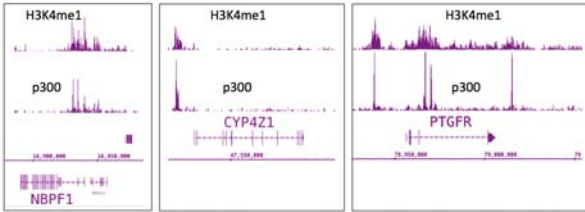
ICC/IF: 0.5 µg/ml dilution

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

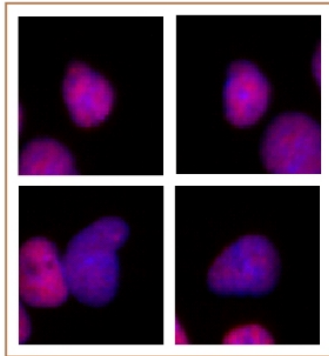
**p300 antibody (mAb) tested by ChIP-Seq.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with chromatin from 4.5 million LNCaP cells and 5 µl of p300 antibody. ChIP DNA was sequenced on the Illumina HiSeq and 15 million sequence tags were mapped to identify p300 binding sites. p300 along with H3K4me1 are markers of active enhancer elements and are therefore expected to co-localize. A sampling of the p300 ChIP-Seq data shows the expected co-localization of p300 and H3K4me1.



**p300 antibody (mAb) tested by ChIP.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with chromatin from 4.5 million LNCaP cells and 4 µg of p300 antibody. ChIP DNA was used in qPCR with the negative control primer pair or gene-specific primer pairs as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.



**p300 antibody (mAb) tested by immunofluorescence.**

Formaldehyde fixed HeLa cells stained with p300 antibody at a 0.5 µg/ml dilution.