

p300 antibody (mAb)

Catalog Nos: 61401, 61903, 61402

RRID: AB_2716754

Clone: NM11

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

Reactivity: Human

Quantities: 100 µg, 50 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Isotype: IgG1

Concentration: 0.1 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.

NOTE: The concentration of this antibody is 0.1 µg/µl.

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

Buffer: Purified IgG in PBS with 50% glycerol.

Application Notes:

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.