

p300 antibody (rAb), 100 µg

Catalog No: 91675**Clone:** 230369D12**Application(s):** IP**Reactivity:** Human**Quantity:** 100 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 264 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: This information is proprietary to Active Motif and/or its suppliers.

Buffer: Purified IgG in PBS, pH 7.3, with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

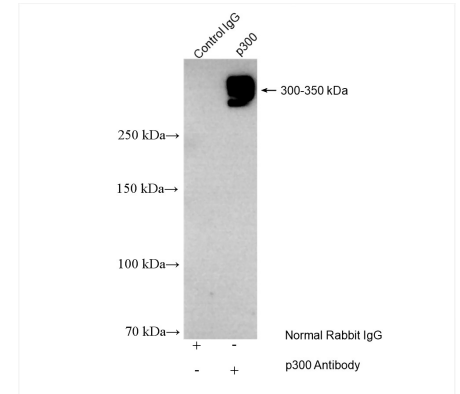
Application Notes:

Applications Validated by Active Motif:

IP: 4 µg per IP

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 (pAb) tested by Immunoprecipitation.**

4 µg of p300 antibody was used to immunoprecipitate p300 from 1050 µg of HeLa nuclear cell extract (lane 2). 10 µl of IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the p300 antibody at a dilution of 1:1000.