

p300 Polyclonal Antibody

(Catalog #A-4020)

Background

p300, also known as EP300 or E1A binding protein p300, is a universally expressed, reasonably stable nuclear phosphorprotein that is closely related to CBP (or CREB). CBP and p300 are identified as transcriptional coactivators that can directly interact with several different DNA-binding transcription factors including nuclear hormone receptors, CREB, c-Fos, C-Jun/v-Jun, c-Myb/v-Myb, TFIIB and MyoD. These two proteins have been known to exhibit histone acetyltransferase (HAT) activity and can acetylate all four core histones in the nucleosome. It is believed that both p300 and CBP may be directly involved in activating chromatin for transcription.

Description

p300 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid in PBS containing 50% glycerol, 0.5% BSA* and 0.02% sodium azide.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q09472

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from Human p300 around the non-phosphorylation site of S89

Storage

Shipped at 4°C. Aliquot store at -20°C. Avoid repeated freeze / thaw cycles.

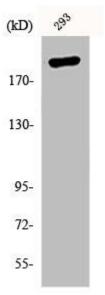
Alternative Names

CREBBP/EP300 inhibitory protein 1 antibody, P-300, Cyclic AMP responsive enhancer binding protein antibody, E1A associated protein p300 antibody, E1A binding protein p300 antibody, EC 2.3.1.48 antibody, EP300 antibody, Histone acetyltransferase p300 antibody, p300 HAT antibody, RB and P300 binding protein EID 1 antibody, Retinoblastoma protein associated protein antibody, RSTS2 antibody

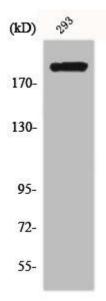
Application

WB, IHC, ELISA; Recommended dilution: WB:1:500-1:2000, IHC:1:100-1:300, ELISA:1:5000

*The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.



Western Blot analysis of 293 cells using p300 Polyclonal Antibody



Western Blot analysis of 293 cells using p300 Polyclonal Antibody