

HP1-alpha Polyclonal Antibody

(Catalog #A-4015)

Background

HP1 α is an isoform of the heterochromatin protein 1 (HP1) which plays a role in chromatin remodeling and transcriptional silencing. The HP1 chromo domain selectively recognizes methylated H3-K9, where this methylation is carried out by histone methyltransferase SUV39H1. SUV39H1-HP1 methylation is important for the correct localization of HP1 at heterochromatin and for transcriptional silencing.

Description

HP1-alpha Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P45973

Purification

Affinity purified

Immunogen

Synthesized peptide derived from the Internal region of human HP1 α .

Storage

Store at -20°C or -80°C. Avoid repeated freeze.

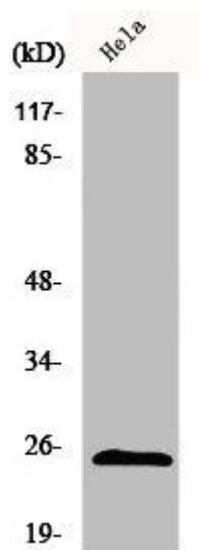
Alternative Names

CBX5; HP1A; Chromobox protein homolog 5; Antigen p25; Heterochromatin protein 1 homolog alpha; HP1 alpha

Application

WB, IHC, IF, ELISA; Recommended dilution: WB:1/500-1/2000, IHC:1/100-1/300, IF:1/200-1/1000, ELISA:1/20000

**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 A549 cells using HP1-alpha Polyclonal Antibody