
Phospho-CDK4-T172 Polyclonal Antibody

(Catalog # A70915)

Background

The protein encoded by this gene is a member of the Ser/Thr protein kinase family. This protein is highly similar to the gene products of *S. cerevisiae* cdc28 and *S. pombe* cdc2. It is a catalytic subunit of the protein kinase complex that is important for cell cycle G1 phase progression. The activity of this kinase is restricted to the G1-S phase, which is controlled by the regulatory subunits D-type cyclins and CDK inhibitor p16(INK4a). This kinase was shown to be responsible for the phosphorylation of retinoblastoma gene product (Rb). Mutations in this gene as well as in its related proteins including D-type cyclins, p16(INK4a) and Rb were all found to be associated with tumorigenesis of a variety of cancers. Multiple polyadenylation sites of this gene have been reported.

Description

Phospho-CDK4-T172 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P11802

Purification

Affinity Purification

Immunogen

A synthetic phosphorylated peptide around T172 of human CDK4 (NP_000066.1).

Storage

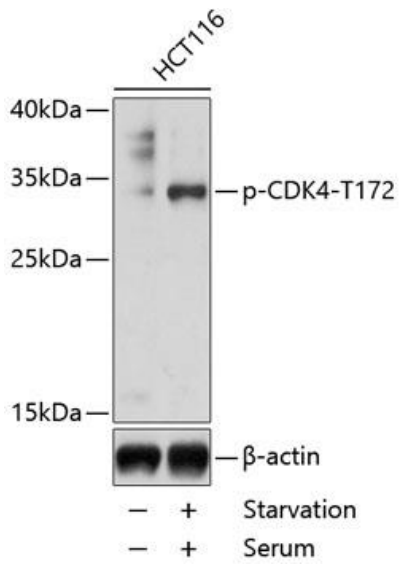
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

CDK4; CMM3; PSK-J3; cyclin-dependent kinase 4

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of extracts of HCT116 cells, using Phospho-CDK4-T172 antibody at 1:1000 dilution. HCT116 cells were treated by 10% FBS for 8 hours after serum-starvation overnight.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% BSA.
Exposure time: 30s.