

# CIPC Polyclonal Antibody

(Catalog # A67674)

## Background

Transcriptional repressor which acts as a negative-feedback regulator of CLOCK-ARNTL/BMAL1 transcriptional activity in the circadian-clock mechanism. Stimulates ARNTL/BMAL1-dependent phosphorylation of CLOCK.

## Description

CIPC Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

## Specificity

Human, Mouse

## Isotype

IgG

## Uniprot ID

Q9C0C6

## Purification

>95%, Protein G purified

## Immunogen

Recombinant Human CLOCK-interacting pacemaker protein (1-399AA)

## Storage

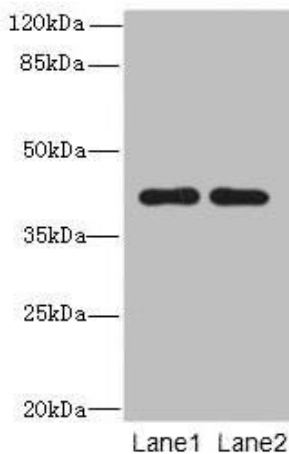
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

## Alternative Names

CLOCK-interacting pacemaker, CLOCK-interacting circadian protein, CIPC, KIAA1737

## Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



## Western blot

All lanes: CIPC Polyclonal Antibody at 6µg/ml

Lane 1: A431 whole cell lysate

Lane 2: NIH/3T3 whole cell lysate

Secondary

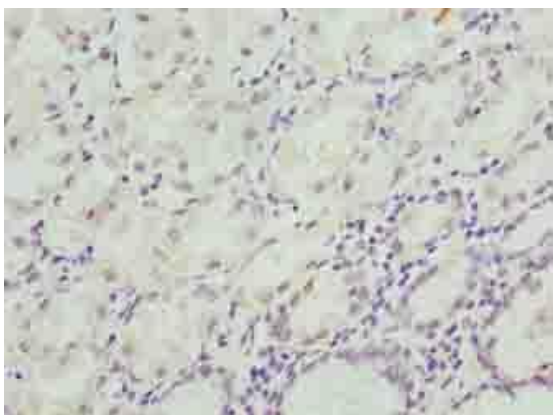
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 43 kDa

Observed band size: 43 kDa



Immunohistochemistry of paraffin-embedded human liver cancer using CIPC Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human stomach tissue using CIPC Polyclonal Antibody at dilution of 1: