
CKAP5 Polyclonal Antibody

(Catalog # A69859)

Background

Binds to the plus end of microtubules and regulates microtubule dynamics and microtubule organization. Acts as processive microtubule polymerase. Promotes cytoplasmic microtubule nucleation and elongation. Plays a major role in organizing spindle poles. In spindle formation protects kinetochore microtubules from depolymerization by KIF2C and has an essential role in centrosomal microtubule assembly independently of KIF2C activity. Contributes to centrosome integrity. Acts as component of the TACC3/ch-TOG/clathrin complex proposed to contribute to stabilization of kinetochore fibers of the mitotic spindle by acting as inter-microtubule bridge. The TACC3/ch-TOG/clathrin complex is required for the maintenance of kinetochore fiber tension. Enhances the strength of NDC80 complex-mediated kinetochore-tip microtubule attachments.

Description

CKAP5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

Q14008

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cytoskeleton-associated protein 5 protein (1802-2019AA)

Storage

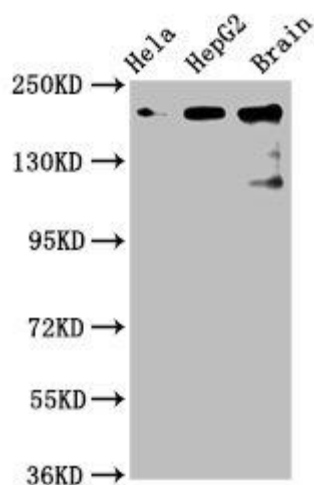
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cytoskeleton-associated protein 5, Colonic and hepatic tumor overexpressed gene protein, Ch-TOG, CKAP5, KIAA0097

Application

ELISA, WB, IHC, IF; Recommended dilution: WB: 1:500-1:5000, IHC: 1:200-1:500, IF: 1:50-1:200



Western Blot

Positive WB detected in: HeLa whole cell lysate, HepG2 whole cell lysate, Rat brain tissue

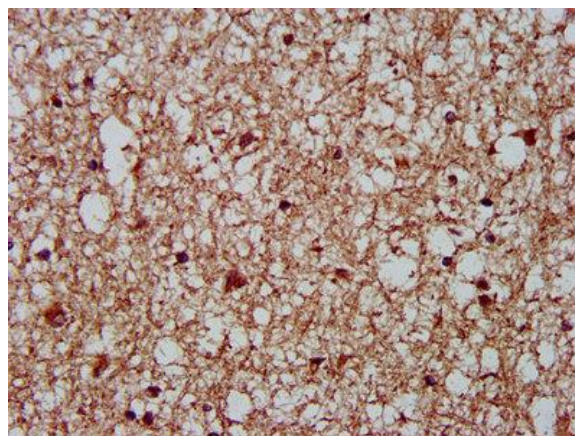
All lanes: CKAP5 Polyclonal Antibody at 6ug/ml

Secondary

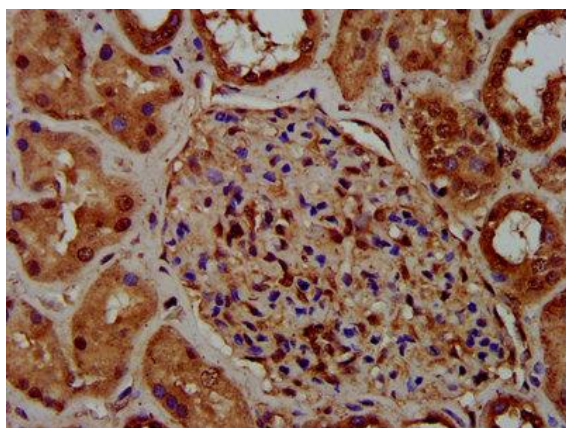
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 226, 219, 227 KDa

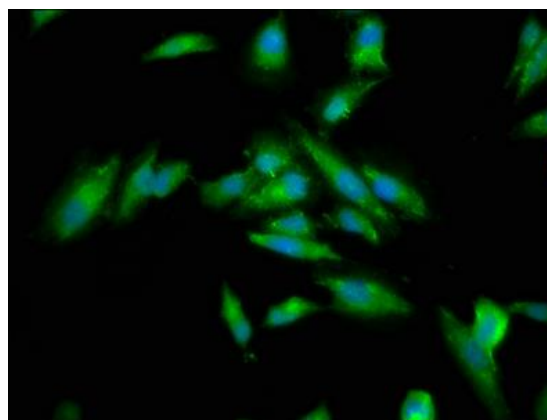
Observed band size: 226 KDa



IHC image of CKAP5 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human brain tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of CKAP5 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HeLa cells with CKAP5 Polyclonal Antibody at 1:133, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).