

CNKSR2 Polyclonal Antibody

(Catalog # A57125)

Background

May function as an adapter protein or regulator of Ras signaling pathways.

Description

CNKSR2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q8WXI2

Purification

Protein G purified

Immunogen

Recombinant Human Connector enhancer of kinase suppressor of ras 2 protein (650-800AA)

Storage

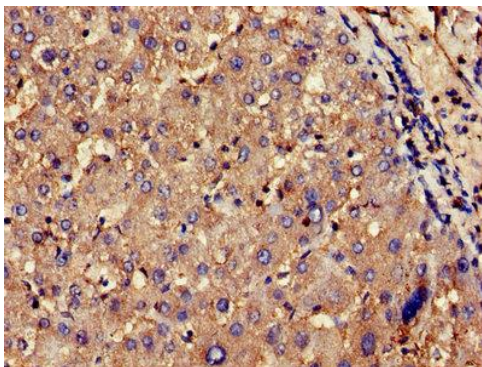
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

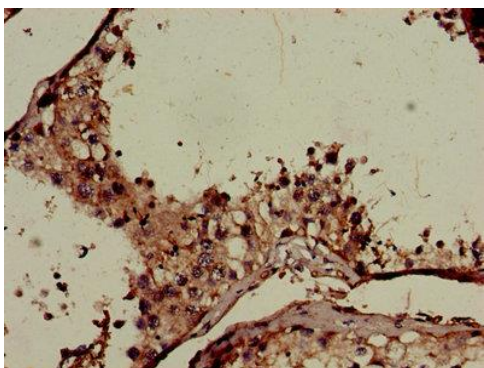
Connector enhancer of KSR 2, CNK homolog protein 2, CNK2, KIAA0902, KSR2, CNKSR2

Application

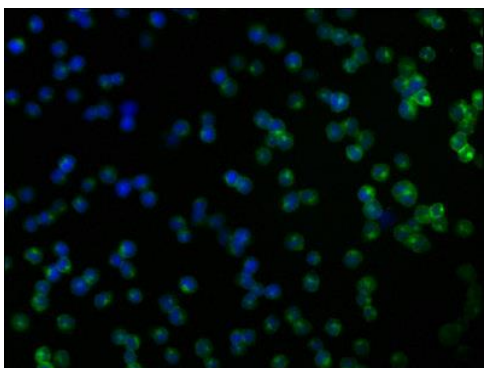
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human liver tissue using CNKSR2 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human testis tissue using CNKSR2 Antibody at dilution of 1:100



Immunofluorescence staining of Jurkat cells with CNKSR2 Antibody at 1:166, 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).