

CDKN2B Polyclonal Antibody

(Catalog # A62250)

Background

Interacts strongly with CDK4 and CDK6. Potent inhibitor. Potential effector of TGF-beta induced cell cycle arrest.

Description

CDKN2B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P42772

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cyclin-dependent kinase 4 inhibitor B protein (1-138AA)

Storage

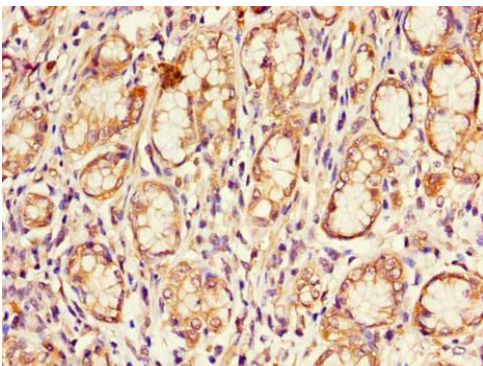
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

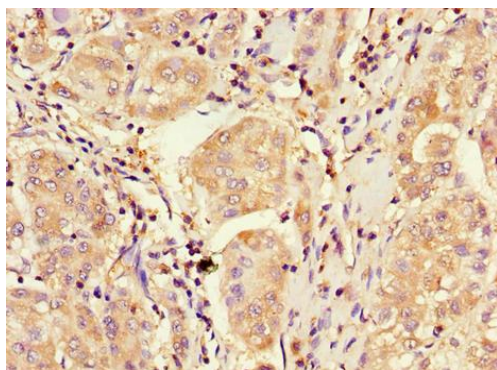
Cyclin-dependent kinase 4 inhibitor B, Multiple tumor suppressor 2, MTS-2, p14-INK4b, p15-INK4b, p15INK4B, CDKN2B, MTS2

Application

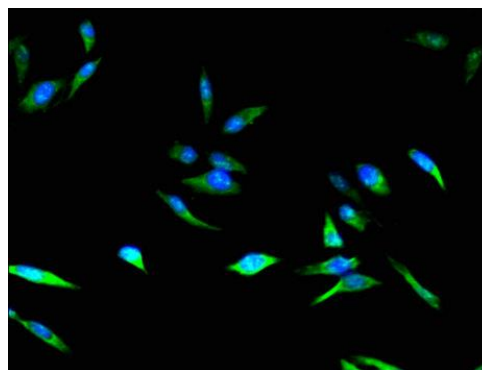
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human gastric tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human liver cancer tissue at dilution of 1:100



Immunofluorescent analysis of HeLa cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)