

CDK2AP1 Polyclonal Antibody

(Catalog #A53256)

Background

specific inhibitor of the cell-cycle kinase CDK2

Description

CDK2AP1 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

O14519

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cyclin-dependent kinase 2-associated protein 1 protein (1-115AA)

Storage

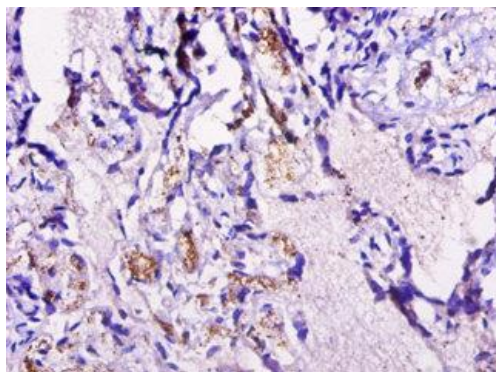
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Deleted in oral cancer 1, Putative oral cancer suppressor, CDK2AP1, CDKAP1, DOC1

Application

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



Immunohistochemistry of paraffin-embedded human placenta at dilution of 1:50