
SYNCRIP Polyclonal Antibody

(Catalog # A51796)

Background

Heterogenous nuclear ribonucleoprotein (hnRNP) implicated in mRNA processing mechanisms. Component of the CRD-mediated complex that promotes MYC mRNA stability. Isoform 1, isoform 2 and isoform 3 are associated in vitro with pre-mRNA, splicing intermediates and mature mRNA protein complexes. Isoform 1 binds to apoB mRNA AU-rich sequences. Isoform 1 is part of the APOB mRNA editosome complex and may modulate the postranscriptional C to U RNA-editing of the APOB mRNA through either by binding to A1CF (APOBEC1 complementation factor), to APOBEC1 or to RNA itself. May be involved in translationally coupled mRNA turnover. Implicated with other RNA-binding proteins in the cytoplasmic deadenylation/translational and decay interplay of the FOS mRNA mediated by the major coding-region determinant of instability (mCRD) domain. Interacts in vitro preferentially with poly(A) and poly(U) RNA sequences. Isoform 3 may be involved in cytoplasmic vesicle-based mRNA transport through interaction with synaptotagmins.

Description

SYNCRIP 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

O60506

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Heterogeneous nuclear ribonucleoprotein Q protein (2-191AA)

Storage

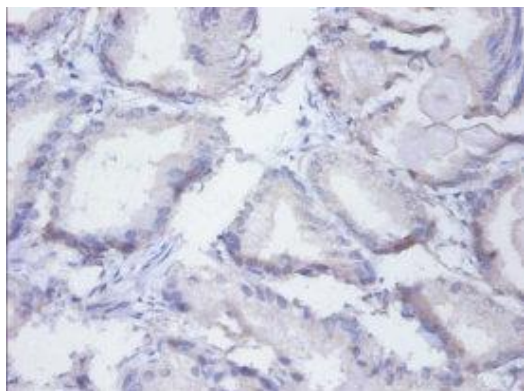
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Application

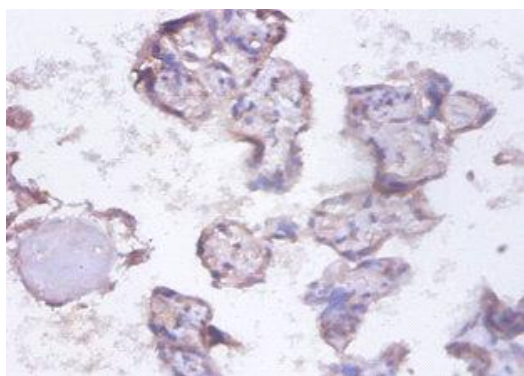
ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



All lanes: SYNCRIP Polyclonal Antibody at 2ug/ml
Lane 1: EC109 whole cell lysate
Lane 2: 293T whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/15000 dilution
Predicted band size: 70, 66, 63, 59, 47 kDa
Observed band size: 69 kDa, 36 kDa



Immunohistochemistry of paraffin-embedded human prostate
at dilution of 1:100



Immunohistochemistry of paraffin-embedded human placenta
at dilution of 1:100