

ROGDI Polyclonal Antibody

(Catalog #A60666)

Background

May act as a positive regulator of cell proliferation.

Description

ROGDI 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

Q9GZN7

Purification

Antigen Affinity Purified

Immunogen

Recombinant Human Protein rogdi homolog protein (14-287AA)

Storage

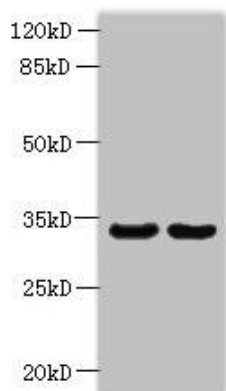
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein rogdi homolog, ROGDI

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200



Western blot

All Lanes: ROGDI Polyclonal Antibody at 1.2ug/ml

Lane 1: Hela whole cell lysate

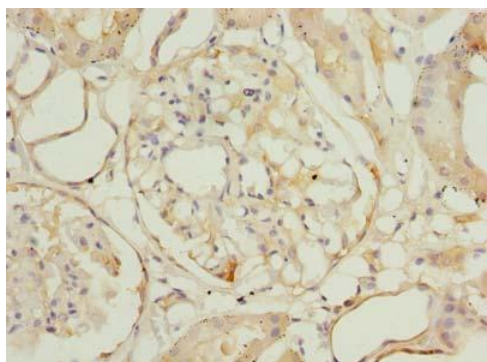
Lane 2: K562 whole cell lysate

Secondary

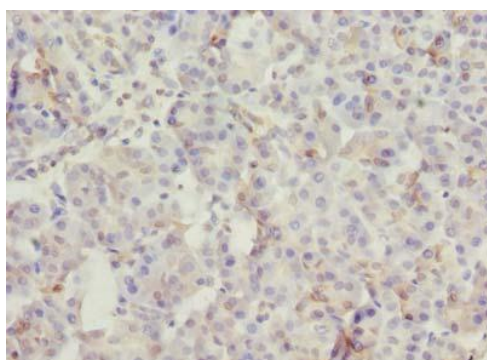
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 33 kDa

Observed band size: 33 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using ROGDI Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human pancreatic tissue using ROGDI Polyclonal Antibody at dilution of 1:100