

C1orf101 Polyclonal Antibody

(Catalog # A65202)

Background

Auxiliary component of the CatSper complex, a complex involved in sperm cell hyperactivation. Sperm cell hyperactivation is needed for sperm motility which is essential late in the preparation of sperm for fertilization.

Description

C1orf101 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q5SY80

Purification

Protein G purified

Immunogen

Recombinant Human Cation channel sperm-associated protein subunit epsilon protein (601-832AA)

Storage

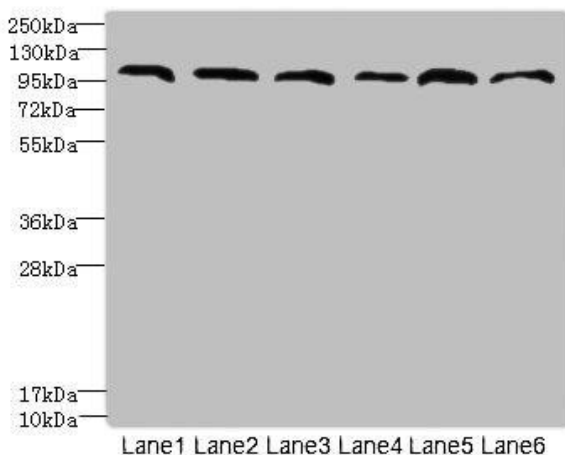
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Uncharacterized protein C1orf101, C1orf101

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All lanes: C1orf101 Polyclonal Antibody at 6ug/ml

Lane 1: U87 whole cell lysate

Lane 2: HepG-2 whole cell lysate

Lane 3: A549 whole cell lysate

Lane 4: Mouse lung tissue

Lane 5: Mouse liver tissue

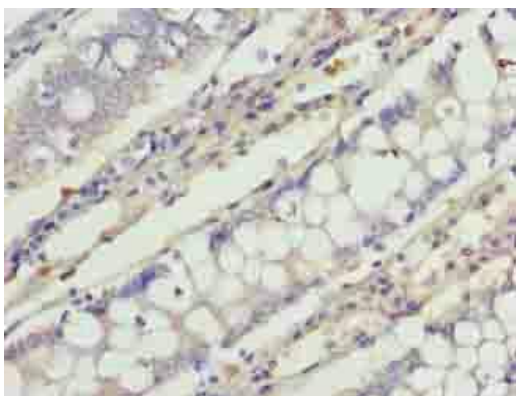
Lane 6: PC-3 whole cell lysate

Secondary

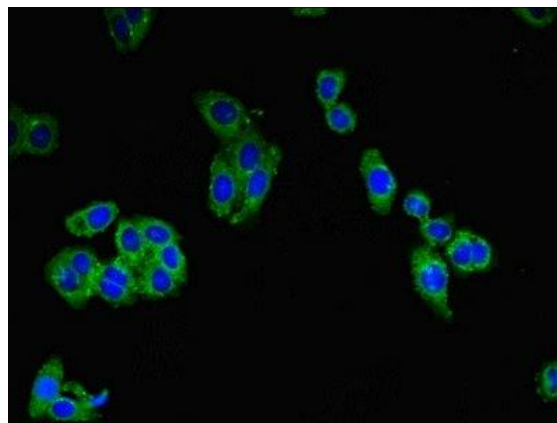
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 110, 96, 93 kDa

Observed band size: 110 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using C1orf101 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using C1orf101 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunohistochemistry of paraffin-embedded human pancreatic cancer using C1orf101 Polyclonal Antibody at dilution of 1:100