

LUZP2 Polyclonal Antibody

(Catalog # A62906)

Description

LUZP2 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

Q86TE4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Leucine zipper protein 2 protein (20-346AA)

Storage

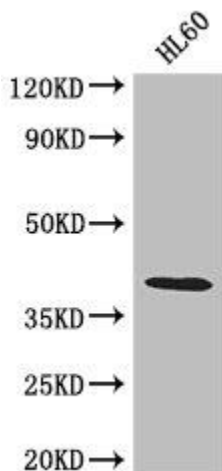
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Leucine zipper protein 2, LUZP2, UNQ2566/PRO6246

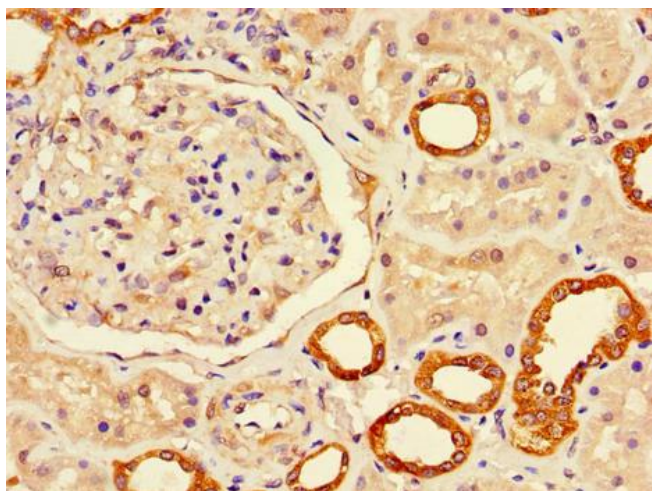
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200

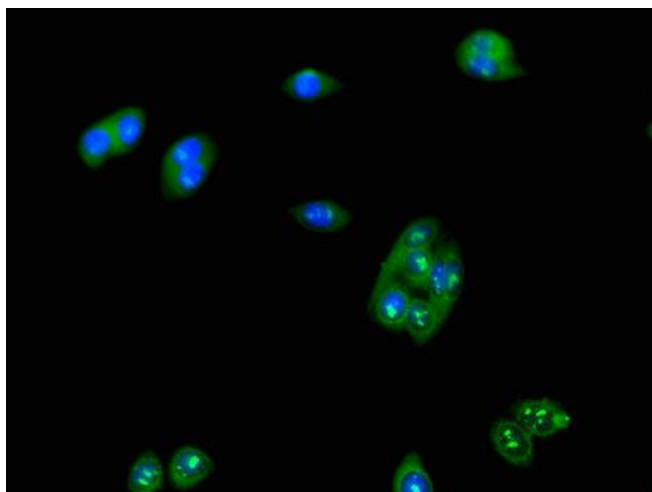


Western Blot

Positive WB detected in: HL6 whole cell lysate
All lanes: LUZP2 Polyclonal Antibody at 2.6ug/ml
Secondary
Goat polyclonal to rabbit IgG at 1/50000 dilution
Predicted band size: 39, 19, 16, 29 kDa
Observed band size: 39 kDa



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



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