

NUP58 Polyclonal Antibody

(Catalog A60106)

Background

Component of the nuclear pore complex, a complex required for the trafficking across the nuclear membrane.

Description

NUP58 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9BVL2

Purification

Protein G purified

Immunogen

Recombinant Human Nucleoporin p58/p45 protein (211-480AA)

Storage

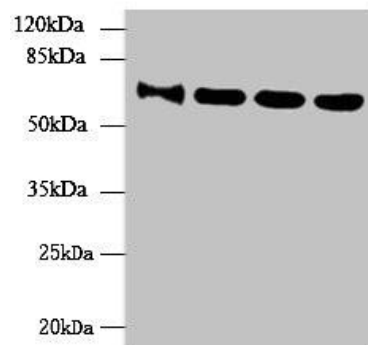
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Nucleoporin p58/p45, NUPL1, KIAA0410, Nucleoporin-like protein 1

Application

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



Western blot

All lanes: NUP58 antibody at 4µg/ml

Lane 1: HepG2 whole cell lysate

Lane 2: Hela whole cell lysate

Lane 3: LO2 whole cell lysate

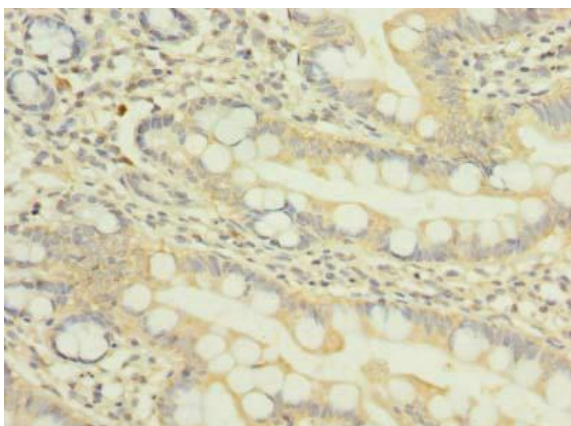
Lane 4: Mouse thymus tissue

Secondary

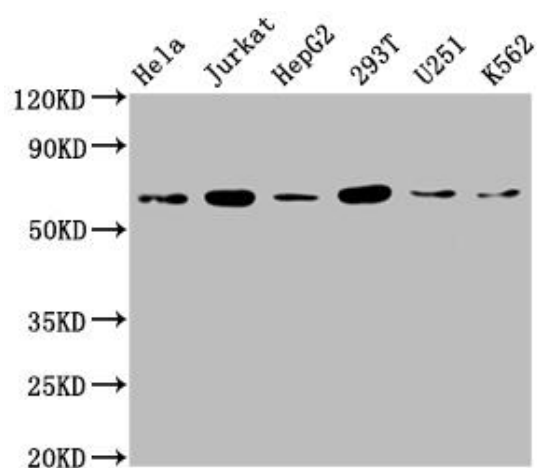
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 61, 51, 60 kDa

Observed band size: 61 kDa



Immunohistochemistry of paraffin-embedded human small intestine tissue using NUP58 Antibody at dilution of 1:100



Western Blot

Positive WB detected in: HeLa whole cell lysate, Jurkat whole cell lysate, HepG2 whole cell lysate, 293T whole cell lysate, U251 whole cell lysate, K562 whole cell lysate

All lanes: NUP58 antibody at 1:2000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 61, 51, 60 kDa

Observed band size: 65 kDa