
NUP85 Polyclonal Antibody

(Catalog # A67538)

Background

Essential component of the nuclear pore complex (NPC) that seems to be required for NPC assembly and maintenance. As part of the NPC Nup107-160 subcomplex plays a role in RNA export and in tethering NUP98/Nup98 and NUP153 to the nucleus. The Nup107-160 complex seems to be required for spindle assembly during mitosis. NUP85 is required for membrane clustering of CCL2-activated CCR2. Seems to be involved in CCR2-mediated chemotaxis of monocytes and may link activated CCR2 to the phosphatidylinositol 3-kinase-Rac-lammellipodium protrusion cascade.

Description

NUP85 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

Q9BW27

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nuclear pore complex protein Nup85 protein (201-500AA)

Storage

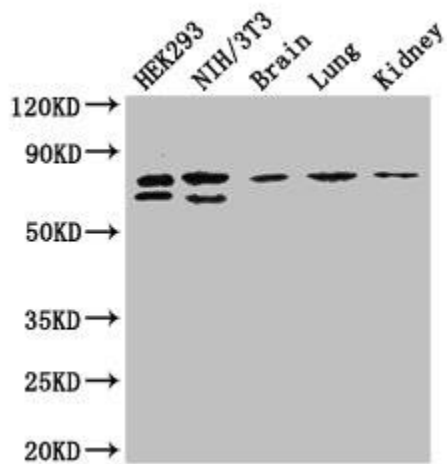
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Nuclear pore complex protein Nup85, 85 kDa nucleoporin, FROUNT, Nucleoporin Nup75, Nucleoporin Nup85, Pericentrin-1, NUP85, NUP75, PCNT1

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: HEK293 whole cell lysate, NIH/3T3 whole cell lysate, Mouse brain tissue, Mouse lung tissue, Mouse kidney tissue

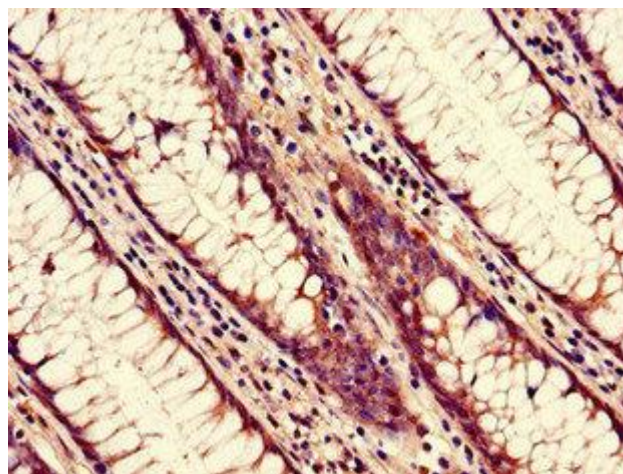
All lanes: NUP85 Polyclonal Antibody at 2.7ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 76, 53, 70 kDa

Observed band size: 76, 70 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using NUP85 Polyclonal Antibody at dilution of 1:100