

## NUP35 Polyclonal Antibody

(Catalog # A66522)

### Background

Functions as a component of the nuclear pore complex (NPC). NPC components, collectively referred to as nucleoporins (NUPs). Can play the role of both NPC structural components and of docking or interaction partners for transiently associated nuclear transport factors. May play a role in the association of MAD1 with the NPC.

### Description

NUP35 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

Q8NFH5

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Nucleoporin NUP53 protein (6-305AA)

### Storage

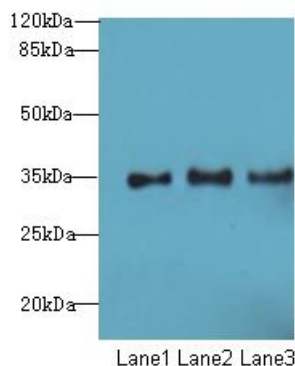
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Nucleoporin NUP53, 35 kDa nucleoporin, Mitotic phosphoprotein 44, MP-44, Nuclear pore complex protein Nup53, Nucleoporin Nup35, NUP35, MP44, NUP53

### Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



#### Western blot

All Lanes: NUP35 Polyclonal Antibody at 12ug/ml

Lane 1: Jurkat whole cell lysate

Lane 2: A431 whole cell lysate

Lane 3: HepG-2 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa