

---

## NUP37 Polyclonal Antibody

(Catalog # A60102)

---

### Background

Component of the Nup107-160 subcomplex of the nuclear pore complex (NPC). The Nup107-160 subcomplex is required for the assembly of a functional NPC. The Nup107-160 subcomplex is also required for normal kinetochore microtubule attachment, mitotic progression and chromosome segregation.

### Description

NUP37 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

Q8NFH4

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Nucleoporin Nup37 protein (1-326AA)

### 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 Nup37, p37, NUP37, Nup107-160 subcomplex subunit Nup37

### Application

ELISA

---