

---

## PELO Polyclonal Antibody, FITC Conjugated

(Catalog #A60232)

---

### Background

Required for normal chromosome segregation during cell division and genomic stability. May function in recognizing stalled ribosomes and triggering endonucleolytic cleavage of the mRNA, a mechanism to release non-functional ribosomes and degrade damaged mRNAs. May have ribonuclease activity (Potential).

### Description

PELO Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9BRX2

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Protein pelota homolog protein (1-385AA)

### 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

Protein pelota homolog (EC:3.1.-.-), PELO, CGI-17

### Application

ELISA; Recommended dilution: ELISA 1:100 - 1:500

---