

---

## ZFP36 Polyclonal Antibody

(Catalog # A68751)

---

### Background

enhances decapping of ARE-containing RNAs, in vitro. Involved in the delivery of target ARE-mRNAs to processing bodies (PBs). In addition to its cytosolic mRNA-decay function, affects nuclear pre-mRNA processing. Negatively regulates nuclear poly(A)-binding protein PABPN1-stimulated polyadenylation activity on ARE-containing pre-mRNA during LPS-stimulated macrophages. Also involved in the regulation of stress granule (SG) and P-body (PB) formation and fusion. Plays a role in the regulation of keratinocyte proliferation, differentiation and apoptosis. Plays a role as a tumor suppressor by inhibiting cell proliferation in breast cancer cells.

### Description

ZFP36 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

P26651

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human mRNA decay activator protein ZFP36 protein (16-103AA)

### Storage

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

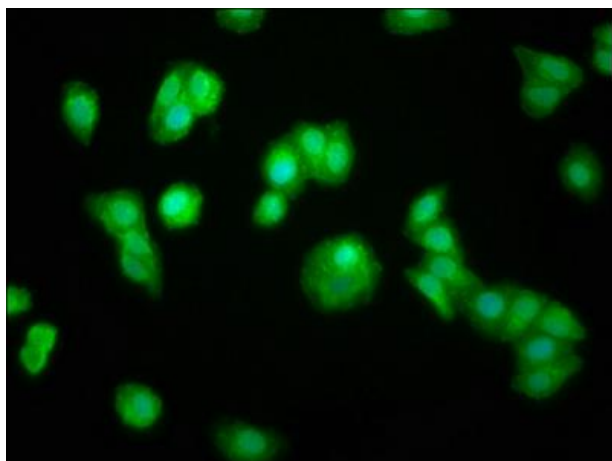
### Alternative Names

mRNA decay activator protein ZFP36, G0/G1 switch regulatory protein 24, Growth factor-inducible nuclear protein NUP475, Tristetraprolin1, ZFP36

### Application

ELISA, IF; Recommended dilution: IF: 1:50-1:200

---



Immunofluorescence staining of HepG2 cells with ZFP36 Polyclonal Antibody at 1:100, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).