

## ZNF98 Polyclonal Antibody

(Catalog # A64048)

### Background

May be involved in transcriptional regulation.

### Description

ZNF98 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

A6NK75

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Zinc finger protein 98 protein (1-135AA)

### Storage

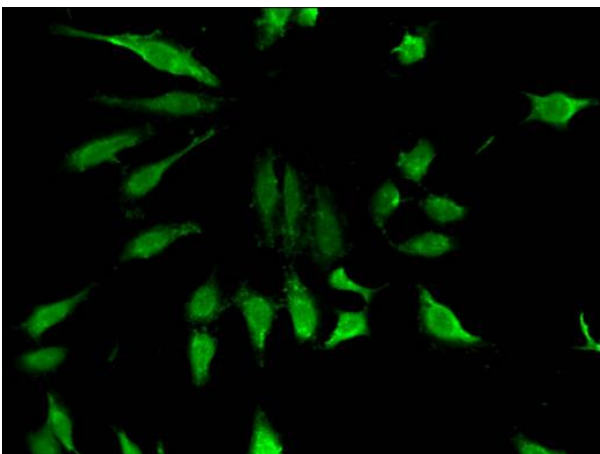
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Zinc finger protein 98, Zinc finger protein 739, Zinc finger protein F7175, ZNF98, ZNF739

### Application

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



Immunofluorescent analysis of HeLa cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)