

## Snrnp35 Polyclonal Antibody

(Catalog # A67810)

### Background

Commitment complex, nucleus, precatalytic spliceosome, U12-type spliceosomal complex, mRNA binding, snRNA binding, mRNA splicing, via spliceosome

### Description

Snrnp35 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q9D384

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Mouse U11/U12 small nuclear ribonucleoprotein 35 kDa protein (1-244AA)

### Storage

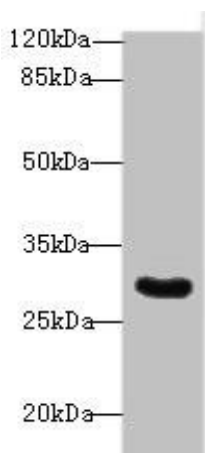
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

U11/U12 small nuclear ribonucleoprotein 35 kDa protein, U11/U12 snRNP 35 kDa protein, U1 snRNP-binding protein homolog, Snrnp35, U1snrnpbp

### Application

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



### Western blot

All lanes: Snrnp35 Polyclonal Antibody at 12ug/ml+Jurkat whole cell lysate

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 29 kDa

Observed band size: 29 kDa