

## Matrix protein Polyclonal Antibody

(Catalog # A54616)

### Background

Plays a major role in assembly and budding of virion. Completely covers the ribonucleoprotein coil and keep it in condensed bullet-shaped form. Inhibits viral transcription and stimulates replication. Plays a major role in early induction of TRAIL-mediated apoptosis in infected neurons

### Description

Matrix protein Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Rabies virus

### Isotype

IgG

### Uniprot ID

P15200

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Rabies virus Matrix protein (1-202AA)

### Storage

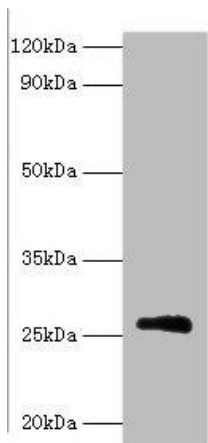
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Phosphoprotein M2 M

### Application

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



### Western blot

All lanes: Matrix protein Polyclonal Antibody at 2ug/ml+recombinant Matrix protein 0.1ug

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 23kDa

Observed band size: 23kDa