

Phosphoprotein Polyclonal Antibody

(Catalog # A54523)

Background

Non catalytic polymerase cofactor and regulatory protein that plays a role in viral transcription and replication. Stabilizes the RNA polymerase L to the N-RNA template and binds the soluble protein N, preventing it from encapsidating non-genomic RNA. Also inhibits host IFN-alpha and IFN-beta signaling by binding and retaining phosphorylated STAT1 in the cytoplasm or by inhibiting the DNA binding of STAT1 in the nucleus. Might be involved, through interaction with host dynein, in intracellular microtubule-dependent virus transport of incoming virus from the synapse toward the cell body

Description

Phosphoprotein 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

P69480

Purification

Protein G purified

Immunogen

Recombinant Rabies virus Phosphoprotein (1-297AA)

Storage

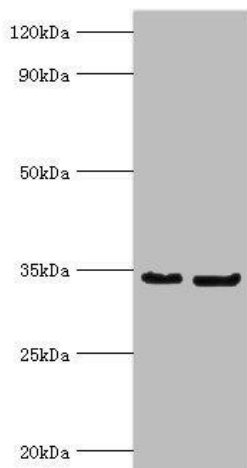
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein M1 P

Application

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



Western blot

All lanes: Phosphoprotein Polyclonal Antibody at 2ug/ml
Lane 1: Rabies virus Phosphoprotein protein at 1ug
Lane 2: Rabies virus Phosphoprotein protein at 0.1ug
Secondary
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 34, 32, 28, 26, 25 kDa
Observed band size: 34 kDa