

Glycoprotein G Polyclonal Antibody

(Catalog # A53464)

Background

Attaches the virus to host cellular receptor, inducing endocytosis of the virion. In the endosome, the acidic pH induces conformational changes in the glycoprotein trimer, which trigger fusion between virus and cell membrane. There is convincing in vitro evidence that the muscular form of the nicotinic acetylcholine receptor (nAChR), the neuronal cell adhesion molecule (NCAM), and the p75 neurotrophin receptor (p75NTR) bind glycoprotein G and thereby facilitate rabies virus entry into cells.

Description

Glycoprotein G 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

P03524

Purification

Protein G purified

Immunogen

Recombinant Rabies virus Glycoprotein protein (20-459AA)

Storage

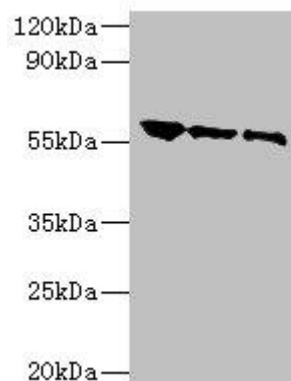
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

G

Application

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



Western blot

All lanes: Glycoprotein G Polyclonal Antibody at 1:500

Lane 1: Rabies virus antigen 1: 2

Lane 2: Rabies virus antigen 1:10

Lane 3: Rabies virus antigen 1:20

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 58kDa

Observed band size: 58kDa