

CD177 Polyclonal Antibody

(Catalog # A66950)

Background

extracellular exosome, plasma membrane, blood coagulation, leukocyte migration

Description

CD177 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

Q8N6Q3

Purification

>95%, Protein G purified

Immunogen

Recombinant Human CD177 antigen protein (22-321AA)

Storage

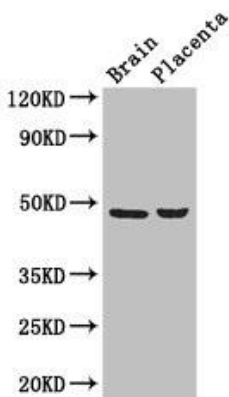
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

CD177 antigen, Human neutrophil alloantigen 2a, HNA-2a, NB1 glycoprotein, NB1 GP, Polycythemia rubra vera protein 1, PRV-1, CD177, CD177, NB1, PRV1, UNQ595/PRO1181

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse brain tissue, Human placenta tissue

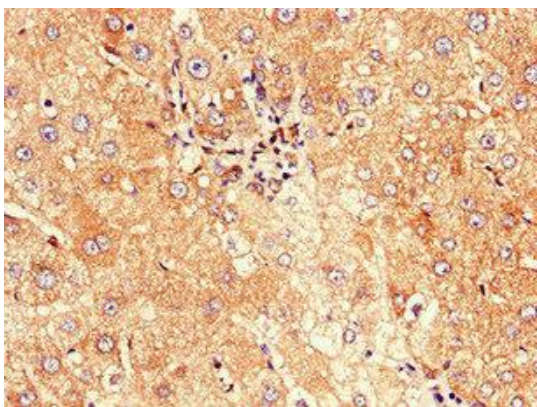
All lanes: CD177 Polyclonal Antibody at 3ug/ml

Secondary

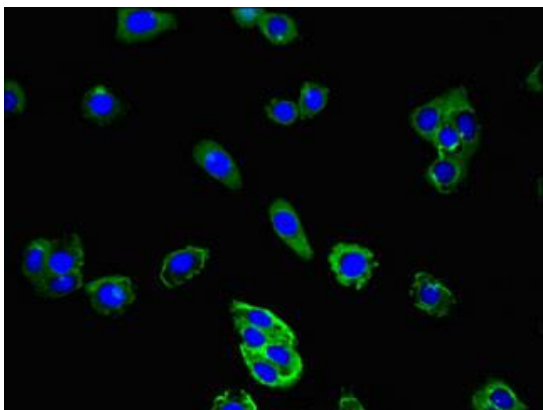
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 47, 28, 16 kDa

Observed band size: 47 kDa



Immunohistochemistry of paraffin-embedded human liver cancer using CD177 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using CD177 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)