
GPA33 Polyclonal Antibody

(Catalog #A51213)

Background

May play a role in cell-cell recognition and signaling.

Description

GPA33 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q99795

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cell surface A33 antigen protein (22-235AA)

Storage

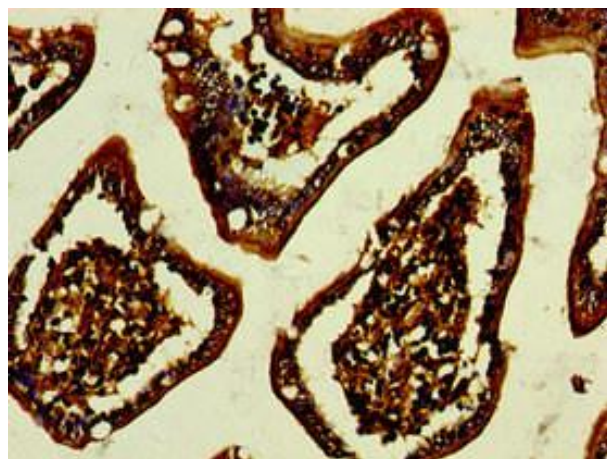
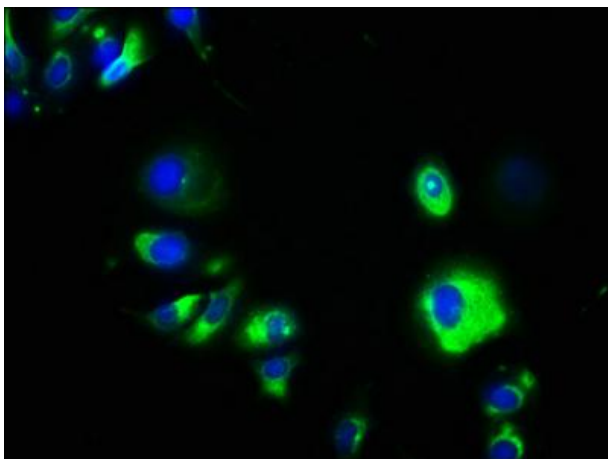
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cell surface A33 antigen, Glycoprotein A33, GPA33

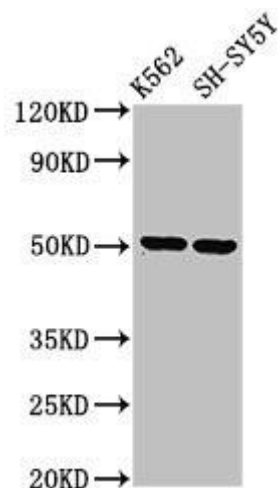
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Immunofluorescence staining of MCF-7 cells with GPA33 Polyclonal Antibody at 1:133, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).

Immunohistochemistry of paraffin-embedded human small intestine tissue using GPA33 Polyclonal Antibody at dilution of 1:100.



Western Blot

Positive WB detected in: K562 whole cell lysate, SH-SY5Y whole cell lysate

All lanes: GPA33 Polyclonal Antibody at 3.2ug/ml
Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 36 KDa

Observed band size: 50 KDa