
GPM6A Polyclonal Antibody

(Catalog # A59282)

Background

Involved in neuronal differentiation, including differentiation and migration of neuronal stem cells. Plays a role in neuronal plasticity and is involved in neurite and filopodia outgrowth, filopodia motility and probably synapse formation. GPM6A-induced filopodia formation involves mitogen-activated protein kinase (MAPK) and Src signaling pathways. May be involved in neuronal NGF-dependent Ca(2+) influx. May be involved in regulation of endocytosis and intracellular trafficking of G-protein-coupled receptors (GPCRs); enhances internalization and recycling of mu-type opioid receptor.

Description

GPM6A 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

P51674

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Neuronal membrane glycoprotein M6-a protein (149-213AA)

Storage

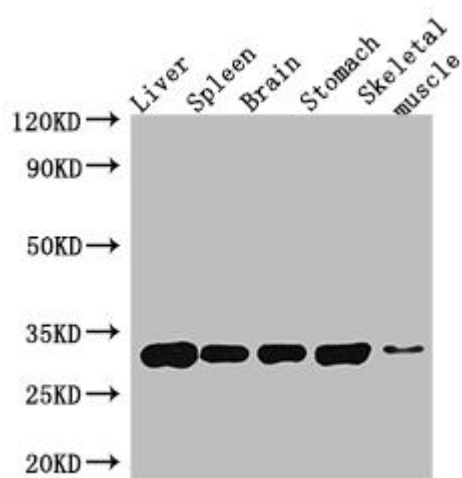
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Neuronal membrane glycoprotein M6-a, M6a, GPM6A, M6A

Application

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



Western Blot

Positive WB detected in: Mouse liver tissue, Mouse spleen tissue, Mouse brain tissue, Mouse stomach tissue, Mouse skeletal muscle tissue

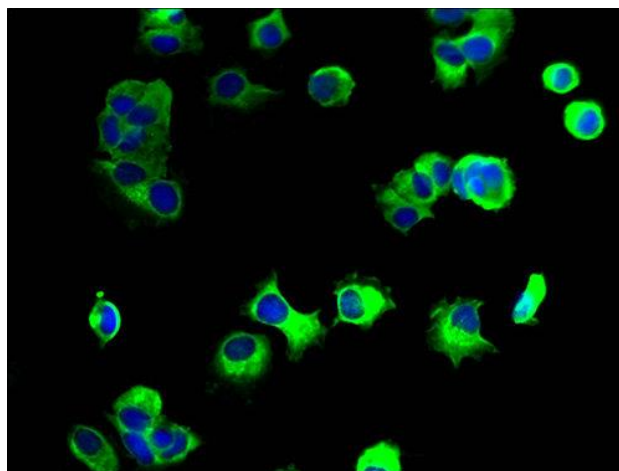
All lanes: GPM6A Polyclonal Antibody at 4ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 32, 31, 30 kDa

Observed band size: 32 kDa



Immunofluorescent analysis of MCF-7 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)