

AMIGO1 Polyclonal Antibody

(Catalog # A65886)

Background

Promotes growth and fasciculation of neurites from cultured hippocampal neurons. May be involved in fasciculation as well as myelination of developing neural axons. May have a role in regeneration as well as neural plasticity in the adult nervous system. May mediate homophilic as well as heterophilic cell-cell interaction and contribute to signal transduction through its intracellular domain. Assembled with KCNB1 modulates the gating characteristics of the delayed rectifier voltage-dependent potassium channel KCNB1.

Description

AMIGO1 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

Q86WK6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Amphoterin-induced protein 1 protein (28-327AA)

Storage

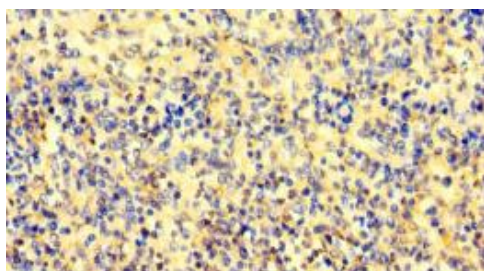
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

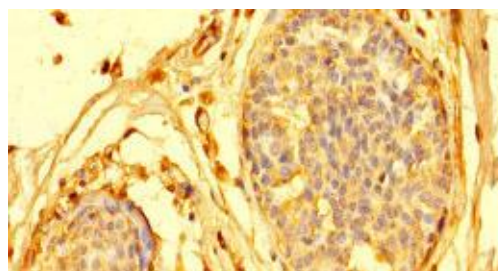
Amphoterin-induced protein 1, AMIGO-1, Alivin-2, AMIGO1

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human spleen tissue using AMIGO1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human breast cancer using AMIGO1 Polyclonal Antibody at dilution of 1:100