

DENND1A Polyclonal Antibody

(Catalog # A58758)

Background

Guanine nucleotide exchange factor (GEF) regulating clathrin-mediated endocytosis through RAB35 activation. Promotes the exchange of GDP to GTP, converting inactive GDP-bound RAB35 into its active GTP-bound form. Regulates clathrin-mediated endocytosis of synaptic vesicles.

Description

DENND1A 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

Q8TEH3

Purification

>95%, Protein G purified

Immunogen

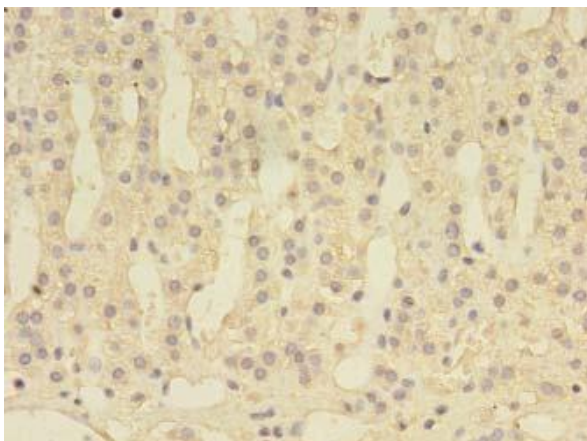
Recombinant Human DENN domain-containing protein 1A protein (1-559AA)

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

DENN domain-containing protein 1A, DENND1A, FAM31A, KIAA1608, Connecdenn 1, Connecdenn, Protein FAM31A



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using DENND1A Polyclonal Antibody at dilution 1:100