
IGSF9 Polyclonal Antibody, FITC Conjugated

(Catalog #A69809)

Background

Functions in dendrite outgrowth and synapse maturation.

Description

IGSF9 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9P2J2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein turtle homolog A protein (132-298AA)

Storage

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

Alternative Names

Protein turtle homolog A, Immunoglobulin superfamily member 9A, IgSF9A, IGSF9, IGSF9A, KIAA1355, NRT1

Application

ELISA; Recommended dilution: ELISA 1:100 - 1:500
