
KIRREL2 Polyclonal Antibody, FITC Conjugated

(Catalog # A65596)

Background

integral component of membrane, plasma membrane, cell adhesion

Description

KIRREL2 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

Q6UWL6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Kin of IRRE-like protein 2 protein (21-219AA)

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

Kin of IRRE-like protein 2, Kin of irregular chiasm-like protein 2, Nephrin-like protein 3, KIRREL2, NEPH3, UNQ5827/PRO19646

Application

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