
KIAA0319L Polyclonal Antibody, FITC Conjugated

(Catalog #A59544)

Background

Possible role in axon guidance through interaction with RTN4R.

Description

KIAA0319L 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

Q8IZA0

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Dyslexia-associated protein KIAA0319-like protein (479-718AA)

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

Dyslexia-associated protein KIAA0319-like protein, KIAA0319L, KIAA1837, PP791

Application

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