
BTNL3 Polyclonal Antibody, FITC Conjugated

(Catalog # A65720)

Description

BTNL3 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

Q6UXE8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Butyrophilin-like protein 3 protein (18-237AA)

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

Butyrophilin-like protein 3, Butyrophilin-like receptor, BTNL3, BTNLR, COLF4100, UNQ744/PRO1472

Application

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