
BTNL9 Polyclonal Antibody, Biotin Conjugated

(Catalog #A65441)

Description

BTNL9 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q6UXG8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Butyrophilin-like protein 9 protein (35-256AA)

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 9, BTNL9, UNQ1900/PRO4346

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
