

BTN2A1 Polyclonal Antibody

(Catalog # A58266)

Background

extracellular exosome, integral component of plasma membrane, lipid metabolic process

Description

BTN2A1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q7KYR7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Butyrophilin subfamily 2 member A1 protein (29-248AA)

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

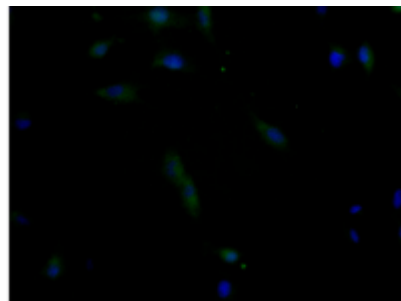
Butyrophilin subfamily 2 member A1, BTN2A1, BT2.1, BTF1

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:300, IF:1:50-1:300



Immunohistochemistry of paraffin-embedded human skeletal muscle tissue using BTN2A1 Polyclonal Antibody at dilution of 1:300



Immunofluorescent analysis of U251 cells using BTN2A1 Polyclonal Antibody at dilution of 1:300 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)