

ATP12A Polyclonal Antibody

(Catalog # A62134)

Background

Catalyzes the hydrolysis of ATP coupled with the exchange of H(+) and K(+) ions across the plasma membrane. Responsible for potassium absorption in various tissues.

Description

ATP12A 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

P54707

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Potassium-transporting ATPase alpha chain 2 protein (171-306AA)

Storage

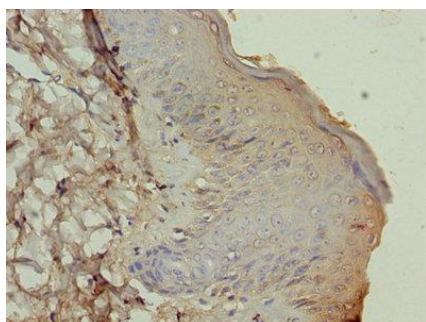
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

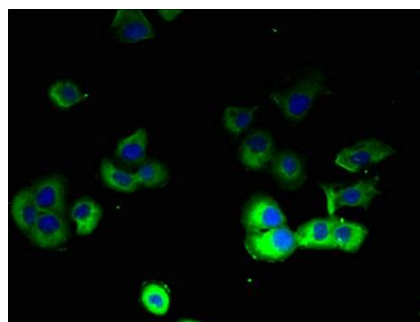
Potassium-transporting ATPase alpha chain 2, Non-gastric H(+)/K(+) ATPase subunit alpha, Proton pump, ATP12A, ATP1AL1

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human skin tissue using ATP12A Antibody at dilution of 1:100



Immunofluorescent analysis of A431 cells using ATP12A Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)