
ATP6V1F Polyclonal Antibody

(Catalog #A52952)

Background

Subunit of the peripheral V1 complex of vacuolar ATPase essential for assembly or catalytic function. V-ATPase is responsible for acidifying a variety of intracellular compartments in eukaryotic cells.

Description

ATP6V1F 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

Q16864

Purification

Protein G purified

Immunogen

Recombinant Human V-type proton ATPase subunit F protein (1-119AA)

Storage

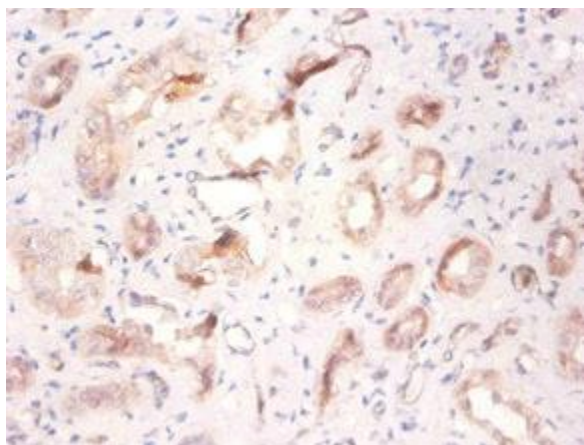
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

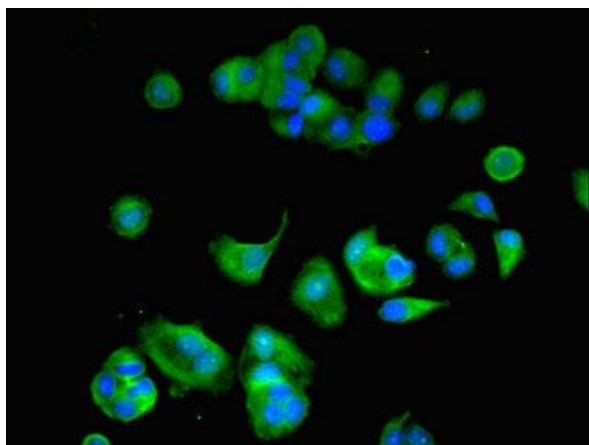
V-ATPase 14 kDa subunit, Vacuolar proton pump subunit F, ATP6V1F, ATP6S14, VATF

Application

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



Immunohistochemistry of paraffin-embedded human kidney tissue using ATP6V1F Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using ATP6V1F Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)