

C9orf72 Polyclonal Antibody

(Catalog # A66710)

Background

May play a role in endosomal trafficking and autophagy.

Description

C9orf72 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

Q96LT7

Purification

Protein G purified

Immunogen

Recombinant Human Guanine nucleotide exchange C9orf72 protein (138-430AA)

Storage

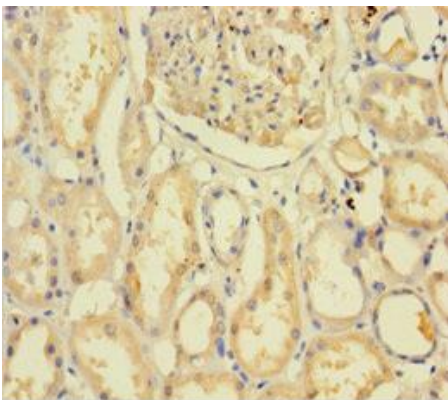
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

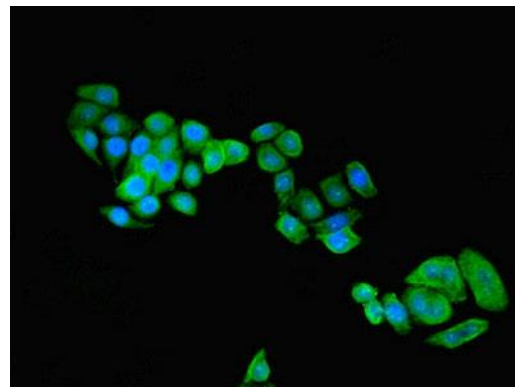
Protein C9orf72, C9orf72

Application

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



Immunohistochemistry of paraffin-embedded human brain tissue using C9orf72 Polyclonal Antibody at dilution 1:100



Immunofluorescent analysis of PC3 cells using C9orf72 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)