

C5orf28 Polyclonal Antibody

(Catalog # A64618)

Description

C5orf28 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

Q0VDI3

Purification

Protein G purified

Immunogen

Recombinant Human Transmembrane protein 267 protein (1-76AA)

Storage

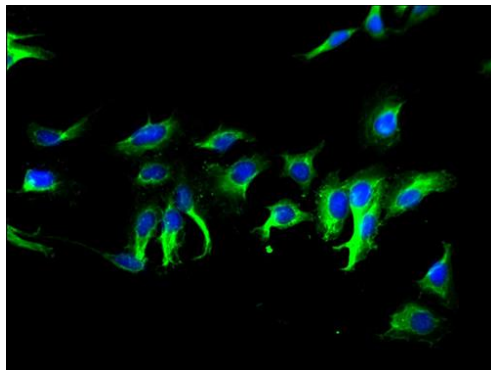
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transmembrane protein C5orf28, C5orf28

Application

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



Immunofluorescent analysis of HeLa cells using C5orf28 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)