

TMEM217 Polyclonal Antibody

(Catalog # A61398)

Description

TMEM217 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

Q8N7C4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane protein 217 protein (150-211AA)

Storage

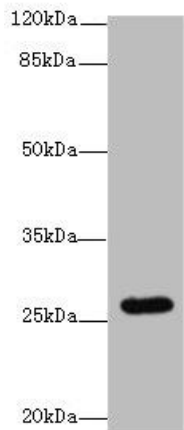
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

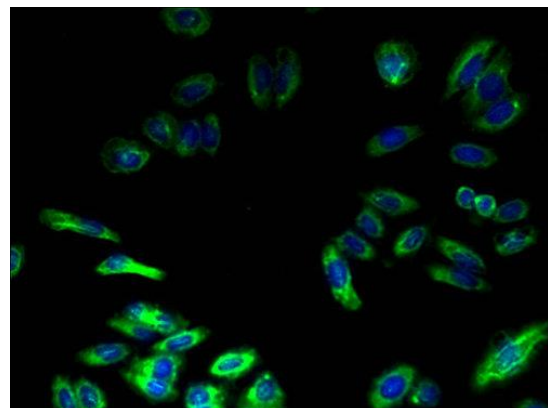
Transmembrane protein 217, TMEM217, C6orf128

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



Western Blot: All lanes-TMEM217 Polyclonal Antibody at 10ug/ml+ Jurkat whole cell lysate
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 27, 25 kDa
Observed band size: 27 kDa



Immunofluorescent analysis of HepG2 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)
