

TMEM222 Polyclonal Antibody

(Catalog # A61402)

Description

TMEM222 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q9H0R3

Purification

Protein G purified

Immunogen

Recombinant Human Transmembrane protein 222 protein (77-164AA)

Storage

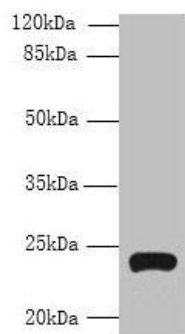
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transmembrane protein 222, TMEM222, C1orf160

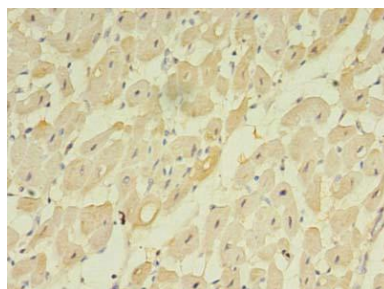
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

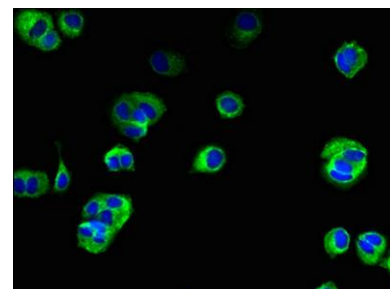


Western blot

All lanes: TMEM222 antibody at 4ug/ml + HepG2: whole cell lysate
Secondary: Goat polyclonal to rabbit IgG at 1/10000 dilution
Predicted band size: 24, 13 kDa
Observed band size: 24 kDa



Immunohistochemistry of paraffin-embedded human heart tissue using TMEM222 Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using TMEM222 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)