

Cxcl5 Polyclonal Antibody

(Catalog # A54093)

Description

Cxcl5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P50228

Purification

Protein G purified

Immunogen

Recombinant Mouse C-X-C motif chemokine 5 protein (46-117AA)

Storage

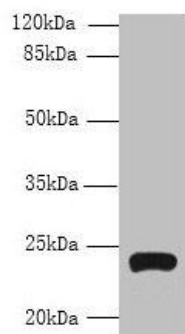
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cytokine LIX Small-inducible cytokine B5

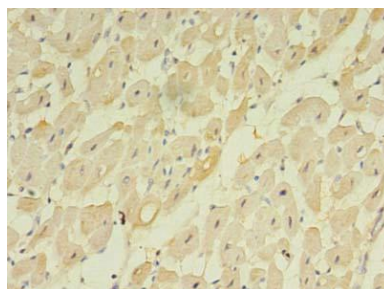
Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000

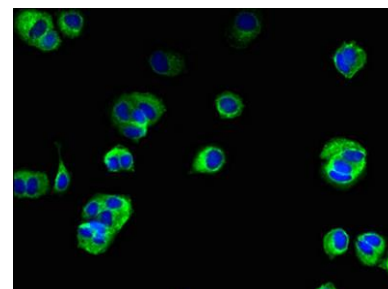


Western blot

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)