

C1QTNF6 Polyclonal Antibody

(Catalog # A67546)

Description

C1QTNF6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9BXI9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Complement C1q tumor necrosis factor-related protein 6 protein (47-278AA)

Storage

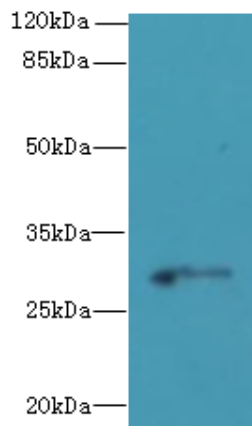
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Complement C1q tumor necrosis factor-related protein 6, C1QTNF6, CTRP6, UNQ581/PRO1151

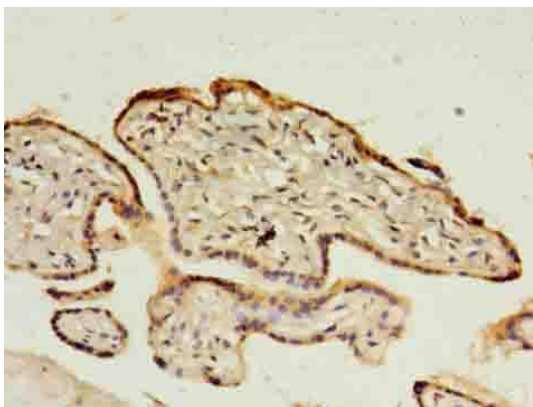
Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200

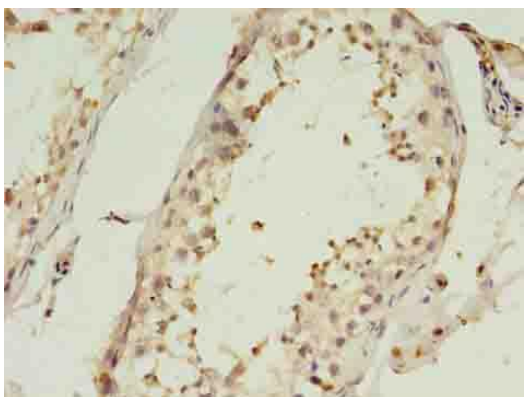


All lanes: C1QTNF6 Polyclonal Antibody at 8ug/ml+Mouse brain tissue
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 29 kDa
Observed band size: 29 kDa

Western blot



Immunohistochemistry of paraffin-embedded human placenta tissue using C1QTNF6 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human testis tissue using C1QTNF6 Polyclonal Antibody at dilution of 1:100