

Plet1 Polyclonal Antibody

(Catalog # A60310)

Background

Modulates leading keratinocyte migration and cellular adhesion to matrix proteins during a wound-healing response and promotes wound repair. May play a role during trichilemmal differentiation of the hair follicle.

Description

Plet1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q8VEN2

Purification

Protein G purified

Immunogen

Recombinant Mouse Placenta-expressed transcript 1 protein (28-218AA)

Storage

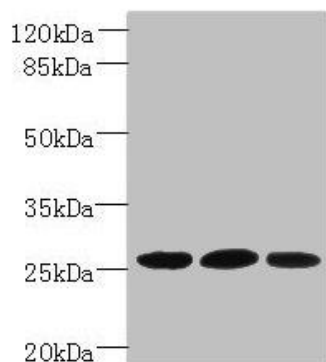
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Placenta-expressed transcript 1 protein, Plet1, Antigen mAgK114

Application

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



Western blot

All lanes: Plet1 antibody at 5ug/ml

Lane 1: L929 whole cell lysate

Lane 2: U251 whole cell lysate

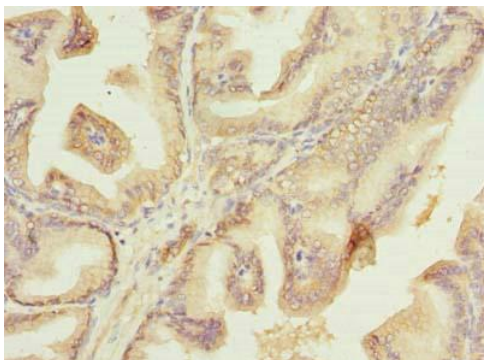
Lane 3: Mouse muscle tissue

Secondary

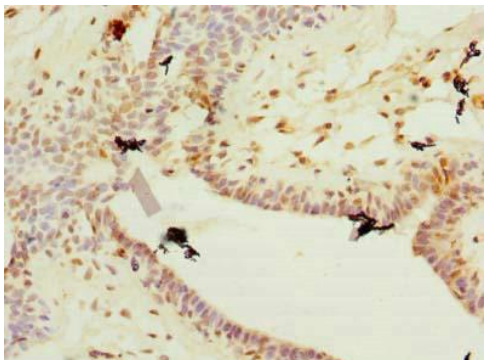
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 26, 21 kDa

Observed band size: 26 kDa



Immunohistochemistry of paraffin-embedded human prostate cancer using Plet1 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human breast cancer using Plet1 Antibody at dilution of 1:100