
MMP3 Polyclonal Antibody

(Catalog # A55223)

Background

Can degrade fibronectin, laminin, gelatins of type I, III, IV, and V; collagens III, IV, X, and IX, and cartilage proteoglycans. Activates procollagenase.

Description

MMP3 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

P08254

Purification

Protein G purified

Immunogen

Recombinant Human Stromelysin-1 protein (100-477AA)

Storage

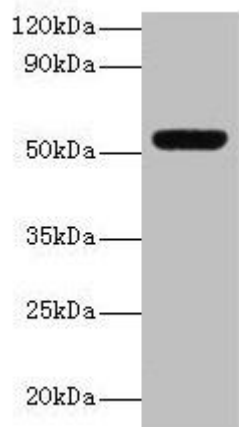
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

SL-1, Matrix metalloproteinase-3, MMP-3, Transin-1, MMP3, STMY1

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200



Western blot

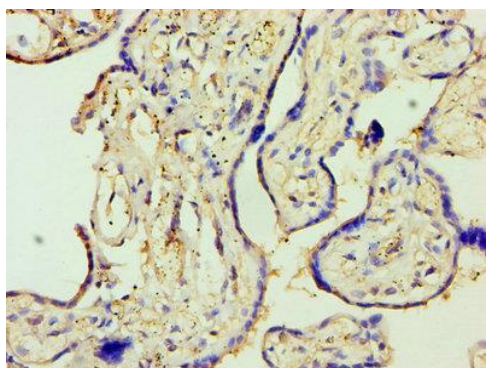
All lanes: MMP3 antibody at 2ug/ml + Mouse liver tissue

Secondary

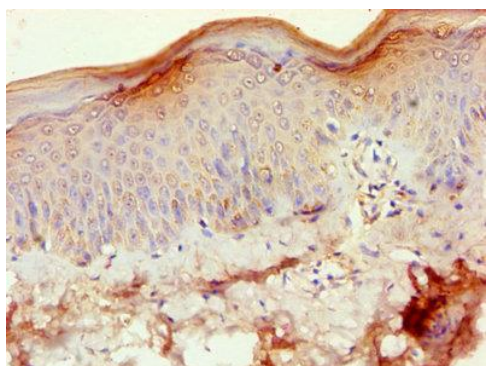
Goat polyclonal to rabbit IgG at 1/15000 dilution

Predicted band size: 54 kDa

Observed band size: 54 kDa



Immunohistochemistry of paraffin-embedded human placenta tissue using MMP3 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human skin tissue using MMP3 Antibody at dilution of 1:100