

MMP10 Polyclonal Antibody

(Catalog #A53300)

Background

Can degrade fibronectin, gelatins of type I, III, IV, and V; weakly collagens III, IV, and V. Activates procollagenase.

Description

MMP10 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Preservative: 0.1mol/L NaCl, 15mmol/L NaN₃, PH7.2.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P09238

Purification

Protein G purified

Immunogen

Recombinant Human Stromelysin-2 protein (99-476AA)

Storage

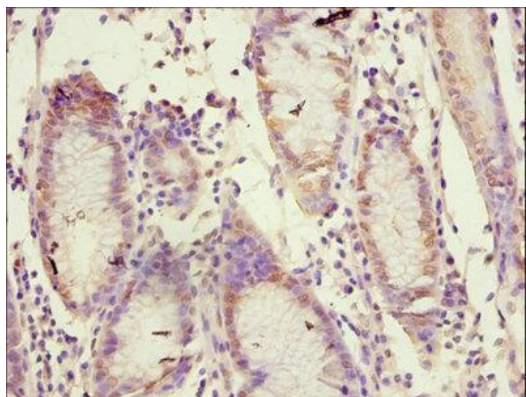
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

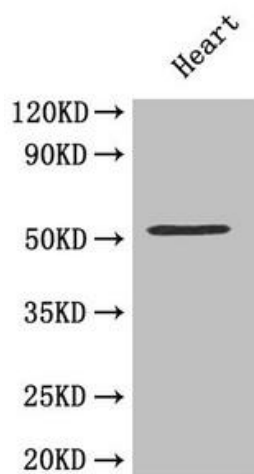
Matrix metalloproteinase-10, Transin-2, MMP10, STMY2

Application

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



Immunohistochemistry of paraffin-embedded human gastric cancer using MMP10 Antibody at dilution of 1:100



Western Blot

Positive WB detected in: Mouse heart tissue

All lanes: MMP10 antibody at 4.5ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 55 kDa

Observed band size: 55 kDa