

---

# HPN Polyclonal Antibody

(Catalog # A69547)

---

## Background

Serine protease that cleaves extracellular substrates, and contributes to the proteolytic processing of growth factors, such as HGF and MST1/HGFL. Plays a role in cell growth and maintenance of cell morphology. Plays a role in the proteolytic processing of ACE2. Mediates the proteolytic cleavage of urinary UMOD that is required for UMOD polymerization.

## Description

HPN Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

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

## Specificity

Human, Mouse, Rat

## Isotype

IgG

## Uniprot ID

P05981

## Purification

>95%, Protein G purified

## Immunogen

Recombinant Human Serine protease hepsin protein (294-413AA)

## Storage

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

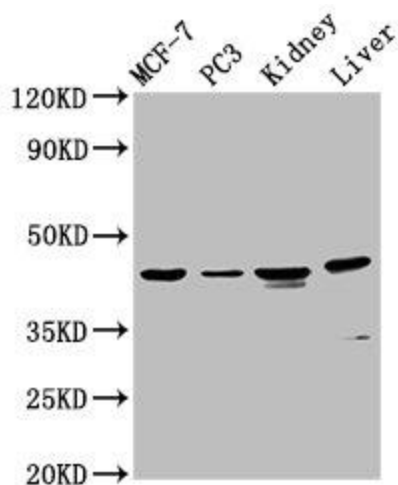
## Alternative Names

Serine protease hepsin, Transmembrane protease serine 1, Serine protease hepsin non-catalytic chain, Serine protease hepsin catalytic chain, HPN, TMPRSS1

## Application

ELISA, WB, IHC, IF; Recommended dilution: WB: 1:500-1:5000, IHC: 1:200-1:500, IF: 1:50-1:200

---



#### Western Blot

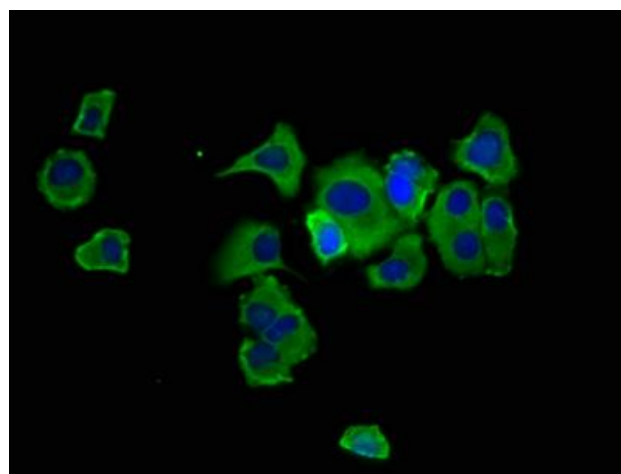
Positive WB detected in: MCF-7 whole cell lysate, PC3 whole cell lysate, Rat kidney tissue, Mouse liver tissue  
All lanes: HPN antibody at 4.5ug/ml

#### Secondary

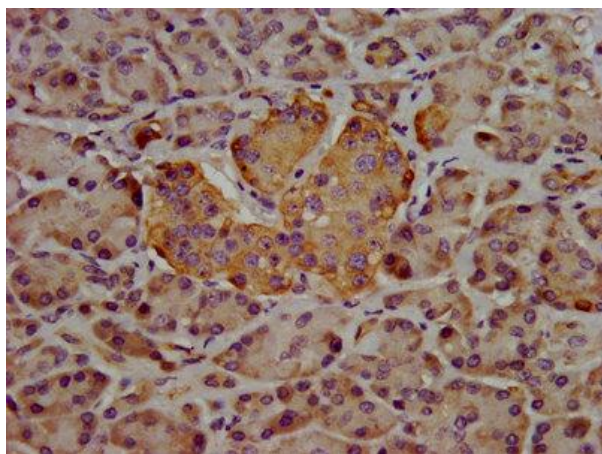
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 46 KDa

Observed band size: 46 KDa



Immunofluorescence staining of MCF-7 cells with HPN Polyclonal Antibody at 1:133, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).



IHC image of HPN Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human pancreatic tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.