
KLK6 Polyclonal Antibody

(Catalog #A66674)

Background

Serine protease which exhibits a preference for Arg over Lys in the substrate P1 position and for Ser or Pro in the P2 position. Shows activity against amyloid precursor protein, myelin basic protein, gelatin, casein and extracellular matrix proteins such as fibronectin, laminin, vitronectin and collagen. Degrades alpha-synuclein and prevents its polymerization, indicating that it may be involved in the pathogenesis of Parkinson disease and other synucleinopathies. May be involved in regulation of axon outgrowth following spinal cord injury. Tumor cells treated with a neutralizing KLK6 antibody migrate less than control cells, suggesting a role in invasion and metastasis.

Description

KLK6 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

Q92876

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Kallikrein-6 protein (22-244AA)

Storage

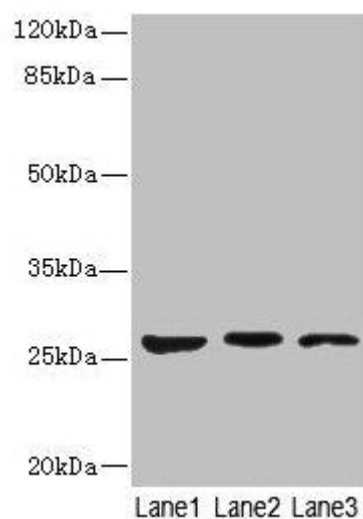
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

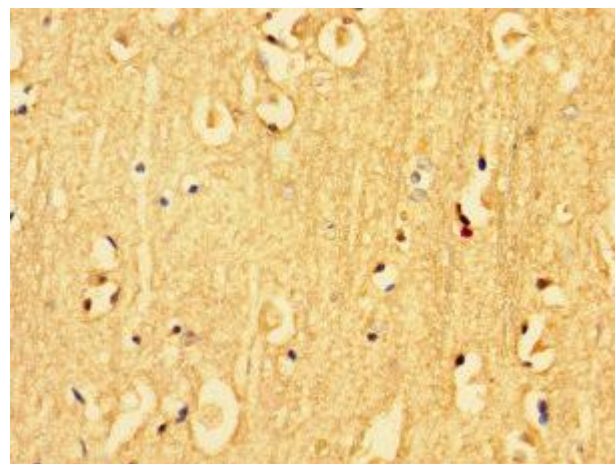
Kallikrein-6, Neurosin, Protease M, SP59, Serine protease 18, Serine protease 9, Zyme, KLK6, PRSS18, PRSS9

Application

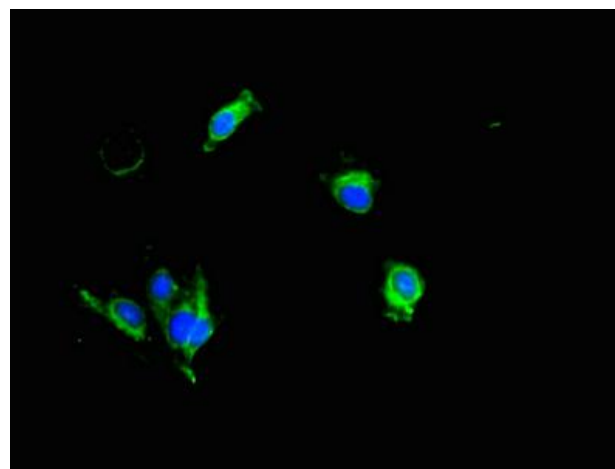
ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



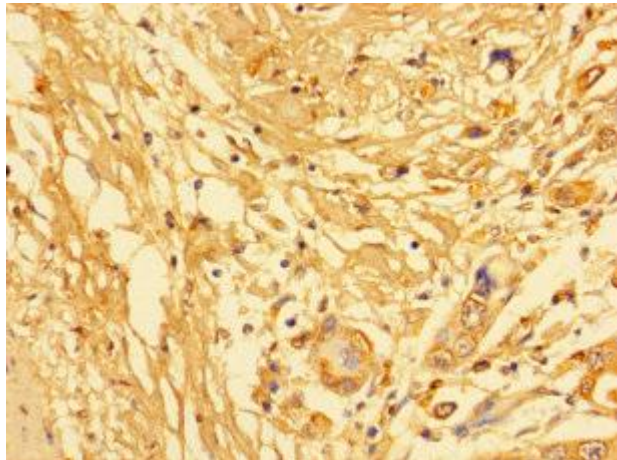
Western blot
 All lanes: KLK6 antibody at 5ug/ml
 Lane 1: Mouse liver tissue
 Lane 2: Mouse brain tissue
 Lane 3: A375 whole cell lysate
 Secondary
 Goat polyclonal to rabbit IgG at 1/10000 dilution
 Predicted band size: 27, 16, 5 kDa
 Observed band size: 27 kDa



Immunohistochemistry of paraffin-embedded human brain tissue using KLK6 Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using KLK6 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human pancreatic cancer using KLK6 Antibody at dilution of 1:100