

## CEP126 Polyclonal Antibody

(Catalog # A55794)

### Background

May participate in cytokinesis.

### Description

CEP126 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9P2H0

### Purification

Protein G purified

### Immunogen

Recombinant Human Centrosomal protein of 126 kDa protein (559-670AA)

### Storage

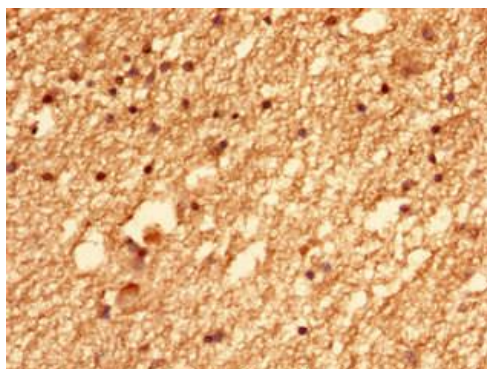
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

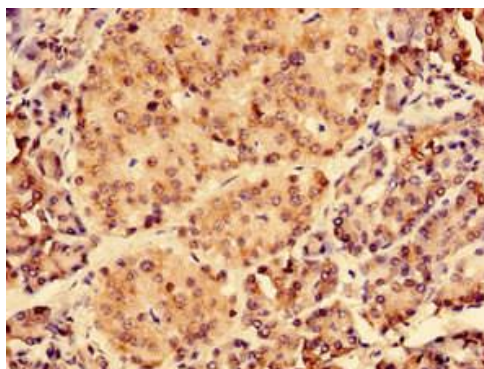
KIAA1377

### Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human brain tissue using CEP126 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human pancreatic tissue using CEP126 Polyclonal Antibody at dilution of 1:100