

KIAA1324L Polyclonal Antibody

(Catalog # A62822)

Description

KIAA1324L 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

A8MWY0

Purification

>95%, Protein G purified

Immunogen

Recombinant Human UPF0577 protein KIAA1324-like protein (227-526AA)

Storage

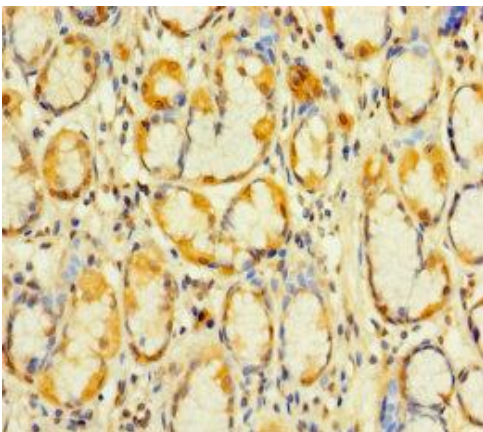
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

UPF0577 protein KIAA1324-like, Estrogen-induced gene 121-like protein, hEIG121L, KIAA1324L, EIG121L

Application

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



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