

LYG1 Polyclonal Antibody

(Catalog # A59698)

Description

LYG1 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

Q8N1E2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Lysozyme g-like protein 1 protein (20-194AA)

Storage

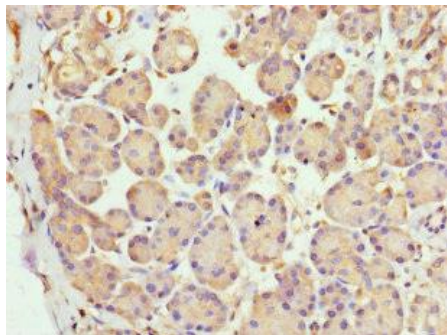
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

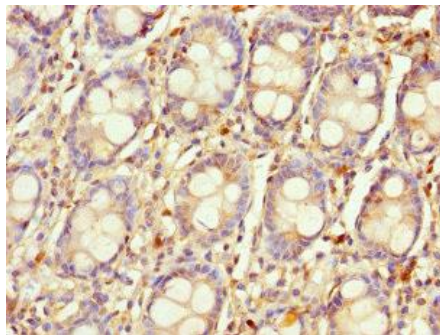
Lysozyme g-like protein 1 (EC:3.2.1.-), LYG1, UNQ1939/PRO4422

Application

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



Immunohistochemistry of paraffin-embedded human pancreatic tissue using LYG1 Polyclonal Antibody at dilution of 1:100. Observed band size: 55kDa, 14kDa



Immunohistochemistry of paraffin-embedded human rectum tissue using LYG1 Polyclonal Antibody at dilution of 1:100