
RER1 Polyclonal Antibody

(Catalog #A60578)

Background

Involved in the retrieval of endoplasmic reticulum membrane proteins from the early Golgi compartment.

Description

RER1 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

O15258

Purification

Protein G purified

Immunogen

Recombinant Human Protein RER1 protein (85-139AA)

Storage

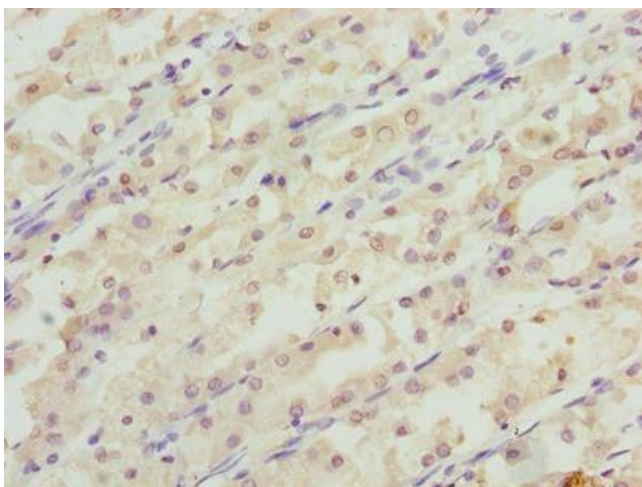
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

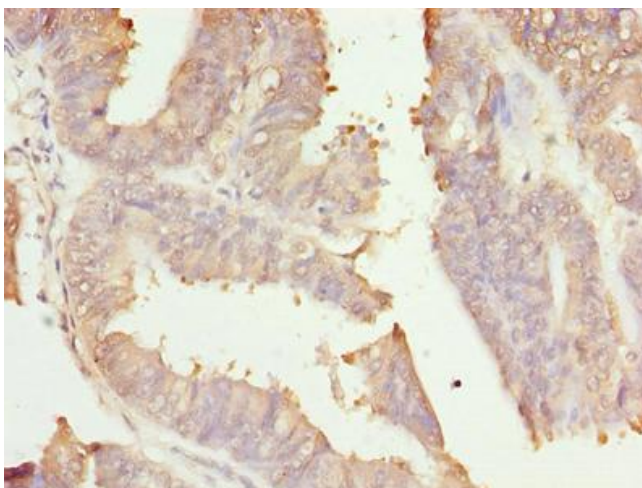
Protein RER1, RER1

Application

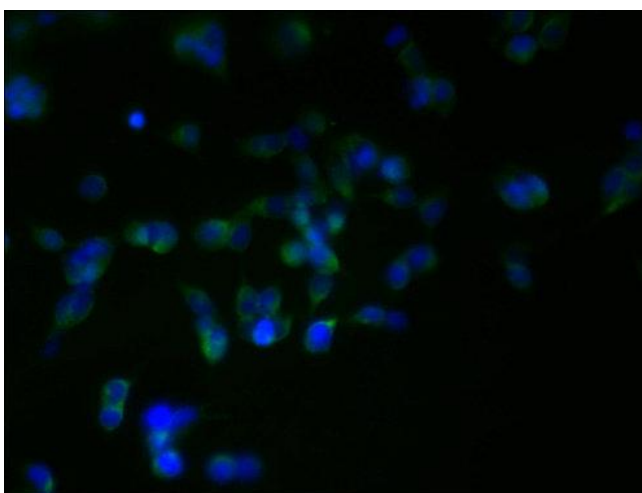
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



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



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



Immunofluorescent analysis of 293T cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)