
RTP4 Polyclonal Antibody

(Catalog # A60742)

Background

Probable chaperone protein which facilitates trafficking and functional cell surface expression of some G-protein coupled receptors (GPCRs). Promotes functional expression of the bitter taste receptor TAS2R16 (PubMed:16720576). Also promotes functional expression of the opioid receptor heterodimer OPRD1-OPRM1.

Description

RTP4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96DX8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Receptor-transporting protein 4 protein (1-224AA)

Storage

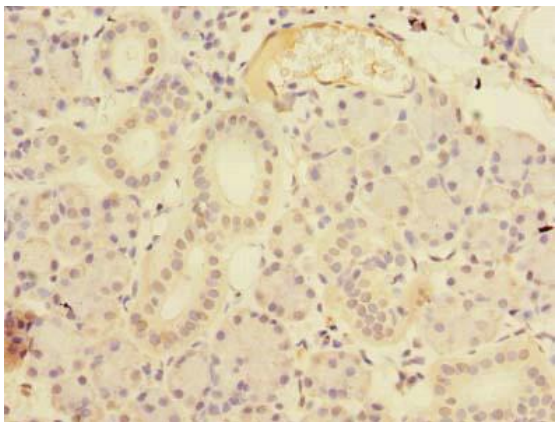
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

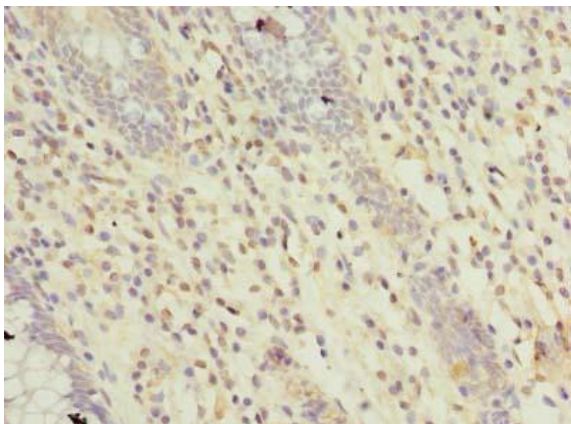
Receptor-transporting protein 4, RTP4, IFRG28, Z3CXXC4, 28 kDa interferon-responsive protein, 3CxxC-type zinc finger protein 4

Application

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



Immunohistochemistry of paraffin-embedded human pancreatic tissue using RTP4 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer using RTP4 Antibody at dilution of 1:100