

GPR183 Polyclonal Antibody

(Catalog # A62698)

Background

Receptor for oxysterol 7- α ,25-dihydroxycholesterol (7- α ,25-OHC) and other related oxysterols. Binding of 7- α ,25-OHC mediates the correct localization of B-cells during humoral immune responses. Promotes activated B-cell localization in the outer follicle and interfollicular regions. Its specific expression during B-cell maturation helps position B-cells appropriately for mounting T-dependent antibody responses. Signals constitutively through G(i)- α , but not through G(s)- α or G(q)- α . Signals constitutively also via MAPK1/3 (ERK1/2).

Description

GPR183 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

P32249

Purification

>95%, Protein G purified

Immunogen

Recombinant Human G-protein coupled receptor 183 protein (1-31AA)

Storage

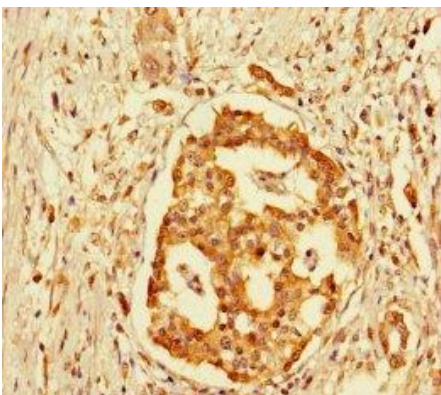
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

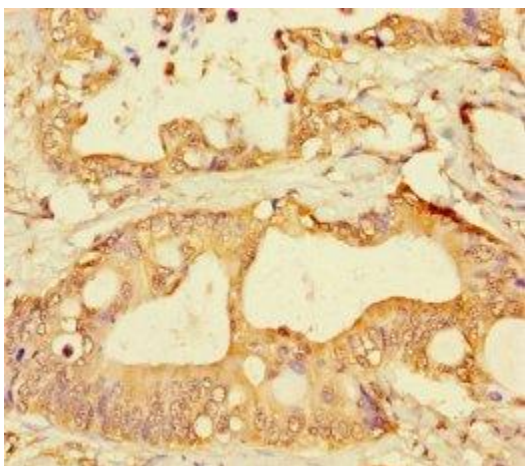
G-protein coupled receptor 183, Epstein-Barr virus-induced G-protein coupled receptor 2, EBI2, EBV-induced G-protein coupled receptor 2, GPR183, EBI2

Application

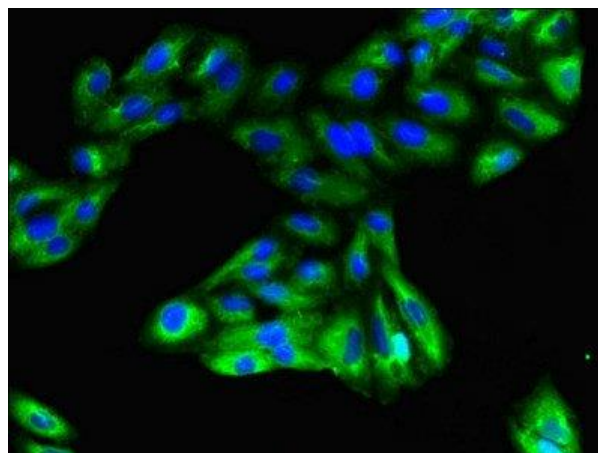
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human pancreatic cancer using GPR183 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer using GPR183 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using GPR183 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)