
GPR84 Polyclonal Antibody

(Catalog #A67298)

Background

Receptor for medium-chain free fatty acid (FFA) with carbon chain lengths of C9 to C14. Capric acid (C10:0), undecanoic acid (C11:0) and lauric acid (C12:0) are the most potent agonists. Not activated by short-chain and long-chain saturated and unsaturated FFAs. Activation by medium-chain free fatty acid is coupled to a pertussis toxin sensitive G(i/o) protein pathway. May have important roles in processes from fatty acid metabolism to regulation of the immune system.

Description

GPR84 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

Q9NQS5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human G-protein coupled receptor 84 protein (202-320AA)

Storage

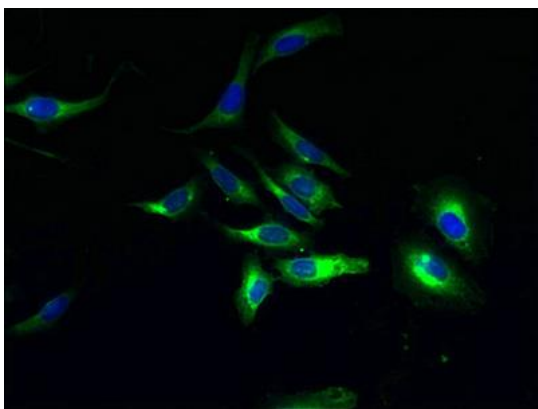
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

G-protein coupled receptor 84, Inflammation-related G-protein coupled receptor EX33, GPR84, EX33

Application

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



Immunofluorescent analysis of HeLa cells using GPR84
Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated
AffiniPure Goat Anti-Rabbit IgG(H+L)