

PRODUCT INFORMATION



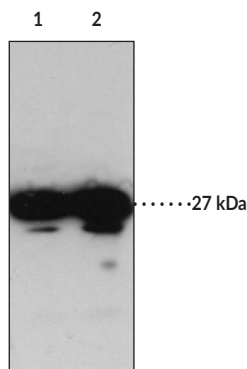
GPR31 (C-Term) Polyclonal Antibody

Item No. 14083

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	G Protein-Coupled Receptor 31, 12(S)-HETE Receptor
Immunogen:	Synthetic peptide from the C-terminal region of human GPR31
Cross Reactivity:	(+) GPR31; (-) GPR41
Species Reactivity:	(+) Human
Uniprot No.:	O00270
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	ELISA, Immunocytochemistry (ICC), and Western blot (WB); the recommended starting dilution for ELISA is 1:500, 1:50 for ICC, and 1:200 for WB. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Image



Lane 1: PC3 cell lysates (40 µg)
Lane 2: PC3 cell lysates (50 µg)

Description

GPR31 is a 319 amino acid G protein-coupled receptor (GPCR) that acts as the receptor for the arachidonic acid metabolite 12-(S)-hydroxy-5,8,10,14-eicosatetraenoic acid (12(S)-HETE) (Item No. 34570). Binding of 12(S)-HETE to GPR31 leads to the activation of ERK1/2 (MAPK3/MAPK1), MEK, and NF-κB pathways. The dysregulation of GPR31 has been implicated in diseases such as arthritis, Alzheimer's disease, and prostate cancer.¹

Reference

1. Guo, Y., Zhang, W., Giroux, C., *et al.* Identification of the orphan G protein-coupled receptor GPR31 as a receptor for 12(S)-hydroxyeicosatetraenoic acid. *J. Biol. Chem.* **286**(39), 33832-33840 (2011).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/13/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM