

PRODUCT INFORMATION



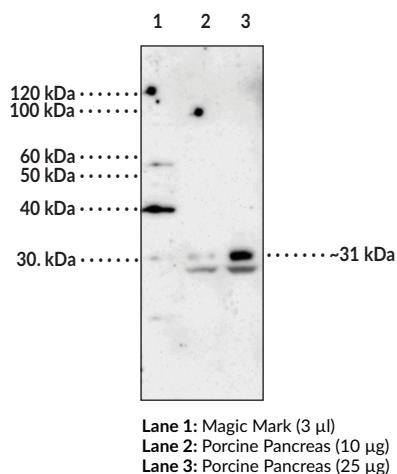
FFAR1 (GPR40) Polyclonal Antibody

Item No. 10007205

Overview and Properties

Contents: This vial contains 500 μ l of peptide affinity-purified polyclonal antibody.
Synonyms: G Protein-Coupled Receptor 40, FFAR1, Free Fatty Acid Receptor
Immunogen: Human GPR40
Species Reactivity: (+) Human, mouse, rat, porcine; other species not tested
Uniprot No.: O14842
Form: Liquid
Storage: -20°C (as supplied)
Stability: \geq 3 years
Storage Buffer: PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host: Rabbit
Applications: Western blot (WB) and immunohistochemistry (IHC) of paraffin-embedded sections; the recommended starting dilution for WB is 1:200 and the recommended starting concentration for IHC is 10 μ g/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Image



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, 10/16/2023

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

PRODUCT INFORMATION



Description

GPR40 is a seven transmembrane G protein-coupled receptor with signalling mediated by medium- and long-chain fatty acids.^{1,2} Found predominantly in the β cells of pancreatic islets, GPR40 has been implicated in the regulation of insulin secretion.^{3,2} Overexpression of GPR40 in murine β cells leads to glucose intolerance suggesting a link between GPR40 expression and diabetes.^{3,6} The receptor has also been identified as a possible mediator of oleate- and linoleate- stimulated proliferation of the human breast cancer cell line MCF-7.^{4,5} GPR40 has a calculated mass of 31.4 kDa based on its amino acid sequence.

References

1. Sawzdargo, M., George, S.R., Nguyen, T., *et al.* A cluster of four novel human G protein-coupled receptor genes occurring in close proximity to CD22 gene on chromosome 19q13.1. *Biochem. Biophys. Res. Commun.* **239**, 543-547 (1997).
2. Itoh, Y., Kawamata, Y., Harada, M., *et al.* Free fatty acids regulate insulin secretion from pancreatic β cells through GPR40. *Nature* **422**, 173-176 (2003).
3. Briscoe, C.P., Peat, A.J., McKeown, S.C., *et al.* Pharmacological regulation of insulin secretion in MIN6 cells through the fatty acid receptor GPR40: Identification of agonist and antagonist small molecules. *Br. J. Pharmacol.* **2006**, 1-10 (2006).
4. Yonezawa, T., Katoh, K., and Obara, Y. Existence of GPR40 functioning in a human breast cancer cell line, MCF-7. *Biochem. Biophys. Res. Commun.* **314**, 805-809 (2004).
5. Hardy, S., St-Onge, G.G., Joly, E., *et al.* Oleate promotes the proliferation of breast cancer cells via the G protein-coupled receptor GPR40. *J. Biol. Chem.* **280**(14), 13285-13291 (2005).
6. Steneberg, P., Rubins, N., Bartoov-Shifman, R., *et al.* The FFA receptor GPR40 links hyperinsulinemia, hepatic steatosis, and impaired glucose homeostasis in mouse. *Cell Metabolism* **1**, 245-258 (2005).

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