

PRODUCT INFORMATION



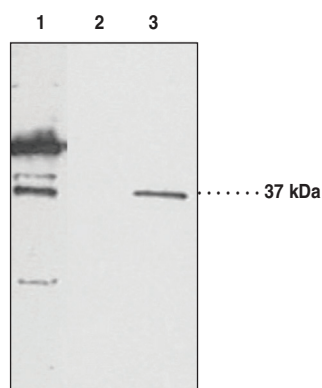
GPR55 Polyclonal Antibody

Item No. 10224

Overview and Properties

Contents:	This vial contains 500 µl of peptide-affinity purified polyclonal antibody.
Synonyms:	G Protein-Coupled Receptor 55, LIP1, Lysophosphatidylinositol Receptor 1
Immunogen:	Synthetic peptide from an internal cytoplasmic region of human GPR55
Species Reactivity:	(+) Human and bovine GPR55
Uniprot No.:	Q9Y2T6
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 and Western blot (WB); the recommended starting dilution is 1:500 and 1:200, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Bovine cornea (5 µg)
Lane 2: HEK293 lysate (13 µg)
Lane 3: GPR55-transfected
HEK293 lysate (13 µg)

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, 11/03/2023

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

GPR55 is an orphan G protein-coupled receptor.¹ GPR55 is expressed in the brain, large dorsal root ganglion neurons, and many peripheral tissues.^{2,3} Endogenous agonists for GPR55 include lysophosphatidylinositol ($EC_{50} = 1.2 \mu\text{M}$ in a β -arrestin-GFP biosensor assay) and the endocannabinoids anandamide (arachidonoyl ethanolamide; Item No. 90050) and 2-arachidonoyl glycerol (Item No. 62160; $EC_{50}s = 18.4$ and 3.5 nM in GTP γ S binding assays).^{1,4} It is also activated by the cannabinoid Δ^9 -tetrahydrocannabinol (Δ^9 -THC; $EC_{50} = 8 \text{ nM}$ in a GTP γ S binding assay).^{1,5} GPR55 is expressed in β cells and pharmacological activation increases glucose-induced insulin release in wild-type mice and, to a lesser extent, in Gpr55 knockout mice.⁶ GPR55 expression is increased in the visceral adipose tissue of obese patients and, to a larger extent, in obese patients with type-2 diabetes.⁷ Activation of GPR55 increases the growth and invasiveness of cancer cells *in vitro*, and its expression in patient-derived tumors is positively correlated with a worse prognosis.⁶ GPR55 activation has also been associated with inhibition of osteoclast formation. Cayman's GPR55 Polyclonal Antibody can be used for flow cytometry, immunofluorescence, and Western blot applications. The antibody recognizes GPR55 at 37 kDa from human and bovine samples. Post-translational modifications such as glycosylation may retard receptor electrophoretic migration such that the protein signal may be detected above 37 kDa.

References

1. Kapur, A., Zhao, P., Sharir, H., *et al.* Atypical responsiveness of the orphan receptor GPR55 to cannabinoid ligands. *J. Biol. Chem.* **284**(43), 29817-29827 (2009).
2. Sawzdargo, M., Nguyenn, T., Lee, D.K., *et al.* Identification and cloning of three novel human G protein-coupled receptor genes GPR52, Ψ GPR53 and GPR55: GPR55 is extensively expressed in human brain. *Brain Res. Mol. Brain Res.* **64**(2), 193-198 (1999).
3. Lauckner, J.E., Jensen, J.B., Chen, H.-Y., *et al.* GPR55 is a cannabinoid receptor that increases intracellular calcium and inhibits M current. *Proc. Natl. Acad. Sci. USA* **105**(7), 2699-2704 (2008).
4. Drmota, T., Greasley, P., and Groblewski, T. Screening assays for cannabinoid-ligand-type modulators of GPR55. (2004).
5. Brown, A.J. Novel cannabinoid receptors. *Br. J. Pharmacol.* **152**, 567-575 (2007).
6. Shore, D.M., and Reggio, P.H. The therapeutic potential of orphan GPCRs, GPR35 and GPR55. *Front. Pharmacol.* **6**, 69, (2015).
7. Moreno-Navarrete, J.M., Catalán, V., Whyte, L., *et al.* The L- α -lysophosphatidylinositol/GPR55 system and its potential role in human obesity. *Diabetes* **61**(2), 281-291 (2012).

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