

GPR55 polyclonal antibody

The cannabinoid receptors CB1 (expressed primarily throughout the central nervous system) and CB2 (expressed in peripheral tissues) are members of the seven-transmembrane domain-containing family of G-protein coupled receptors (GPCRs), and are named for their signaling function in response to cannabinoid drugs such as delta-9-tetrahydrocannabinol (delta-9-THC). Significant evidence suggests that additional cannabinoid receptors may contribute to the many behavioral, vascular, and immunological effects of agonists such as THC and endogenous cannabinoids. Evidence suggests GPR55 may act as a cannabinoid receptor, although its definition as a bona fide "CB3" receptor remains controversial. GPR55 is highly expressed in large-diameter dorsal root ganglion (DRG) neurons, and its activation by cannabinoid compounds (THC, JWH015, and anandamide) has been shown to release calcium from IP3-sensitive ER stores. These findings are opposed by evidence that suggests GPR55 is an intrinsic receptor for lysophosphatidylinositol (LPI).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

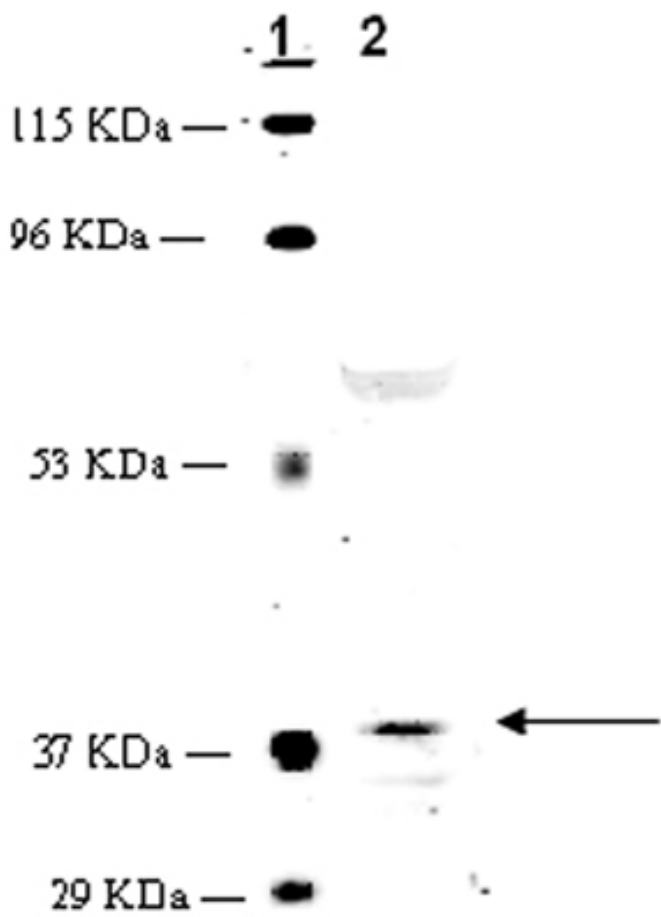
Ordering Information

[Order Online »](#)

ADI-905-900-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW marker (1) and 25 μ g rat kidney extract (2), probed with GPR55/Cannabinoid Receptor, pAb.

Handling & Storage

Handling Keep on ice at all times.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CNR, CBR, GPR55 Cannabinoid Receptor

Application WB

Application Notes Predicted MW of ~34kDa, observed by Western blot ~37kDa. Higher MW species (~60kDa) and intermediates also observed by Western blot which reflect post-translational modification of the receptor.

Formulation Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of rat GPR55.

Purity Detail Peptide affinity purified.

Species Reactivity Mouse, Rat

UniProt ID Q3UJF0 (mouse)

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com