

# 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 »](#)

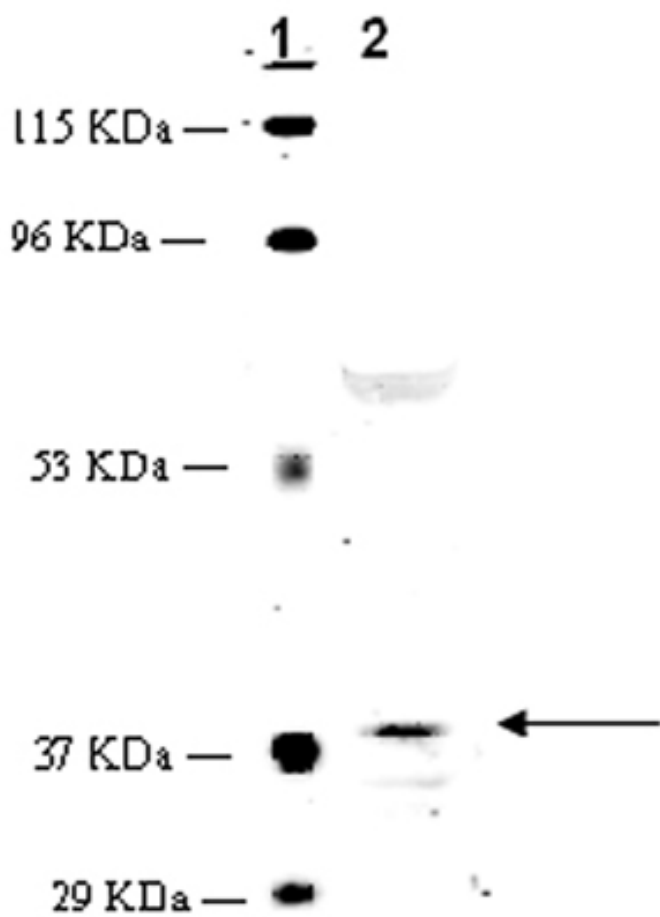
## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW marker (1) and 25µg 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](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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