## **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** 

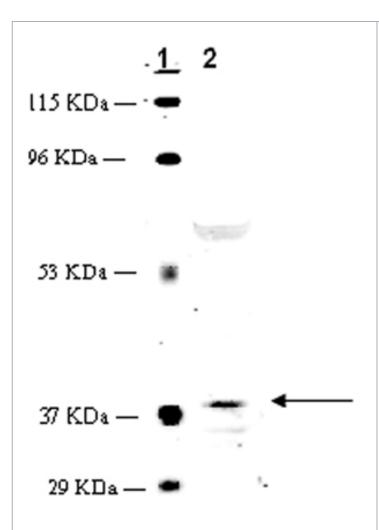
Order Online »

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

•

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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