
MRGPRX2 Polyclonal Antibody

(Catalog # A66606)

Background

Mast cell-specific receptor for basic secretagogues, i. e. cationic amphiphilic drugs, as well as endo- or exogenous peptides, consisting of a basic head group and a hydrophobic core (PubMed:25517090). Recognizes and binds small molecules containing a cyclized tetrahydroisoquinoline (THIQ), such as non-steroidal neuromuscular blocking drugs (NMBDs), including tubocurarine and atracurium. In response to these compounds, mediates pseudo-allergic reactions characterized by histamine release, inflammation and airway contraction. Acts as a receptor for a number of other ligands, including peptides and alkaloids, such as cortistatin-14, proadrenomedullin N-terminal peptides PAMP-12 and, at lower extent, PAMP-20, antibacterial protein LL-37, PMX-53 peptide, beta-defensins, and compleanadine A.

Description

MRGPRX2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96LB1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Mas-related G-protein coupled receptor member X2 protein (286-330AA)

Storage

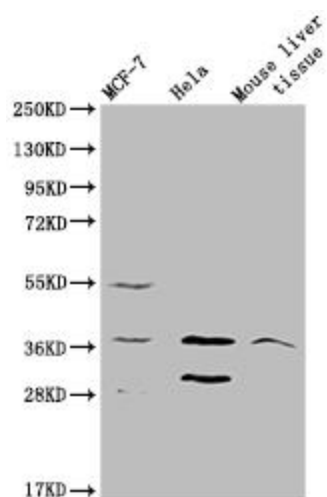
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Mas-related G-protein coupled receptor member X2, MRGPRX2, MRGX2

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



Positive WB detected in: MCF-7 whole cell lysate, HeLa whole cell lysate, Mouse liver tissue

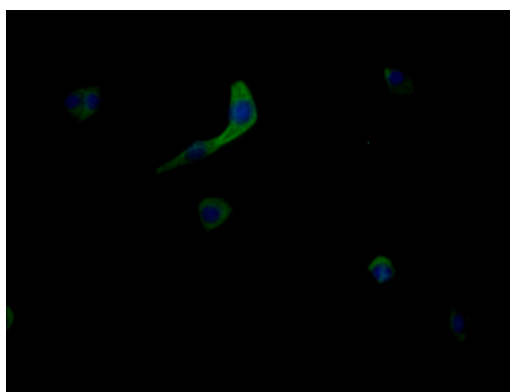
All lanes: MRGPRX2 Polyclonal Antibody at 1:2000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 38 kDa

Observed band size: 38 kDa



Immunofluorescent analysis of MCF-7 cells using MRGPRX2 Polyclonal Antibody at dilution of 1:200 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)