

NMUR1 Polyclonal Antibody

(Catalog #A67838)

Background

Receptor for the neuromedin-U and neuromedin-S neuropeptides.

Description

NMUR1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9HB89

Purification

Protein G purified

Immunogen

Recombinant Human Neuromedin-U receptor 1 protein (360-426AA)

Storage

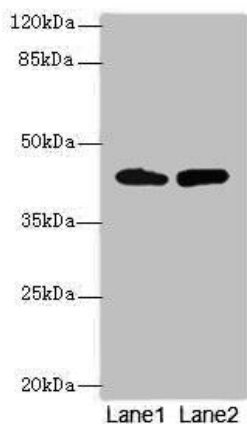
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Neuromedin-U receptor 1, NMU-R1, G-protein coupled receptor 66, G-protein coupled receptor FM-3, NMUR1, GPR66

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

All lanes: NMUR1 antibody at 2ug/ml

Lane 1: Mouse liver tissue

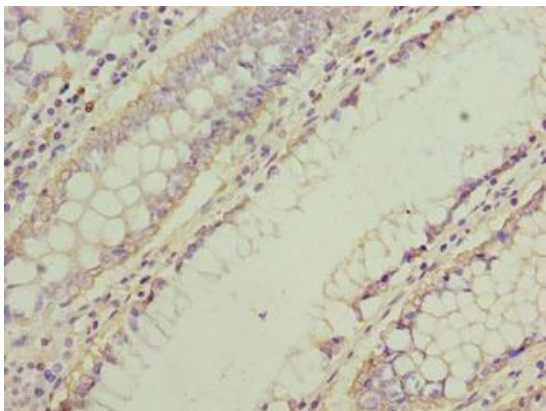
Lane 2: Mouse kidney tissue

Secondary

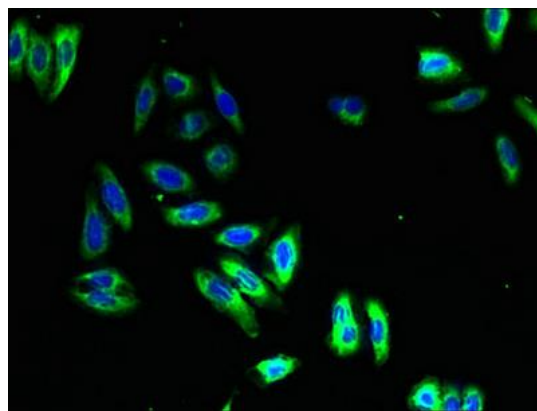
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 47 kDa

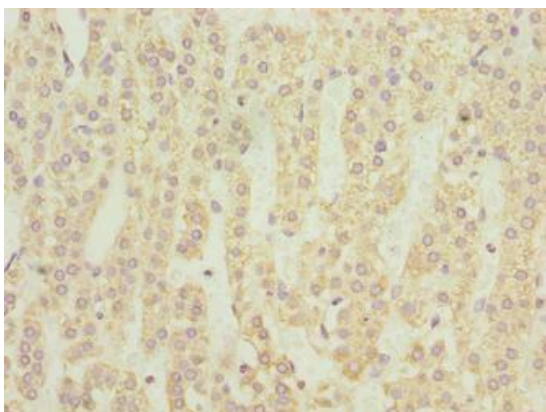
Observed band size: 47 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using NMUR1 Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using NMUR1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using NMUR1 Antibody at dilution of 1:100