

OR2L13 Polyclonal Antibody

(Catalog # A66351)

Background

OR2L13 (Olfactory Receptor Family 2 Subfamily L Member 13) is an odorant receptor protein-coding gene. To initiate a neuronal response that triggers the perception of smell, olfactory receptors interact with odorant molecules in the nose. Olfactory receptor proteins belong to the large family of G-protein-coupled receptors (GPCR) which arise from single coding-exon genes. The olfactory receptor gene family is the largest in the genome. Olfactory receptors share a 7-transmembrane domain structure with many neurotransmitters and hormone receptors. They are responsible for the recognition and G protein-mediated transduction of odorant signals. The nomenclature assigned to the olfactory receptor genes and proteins for this organism is independent of other organisms.

Description

OR2L13 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

Q8N349

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Olfactory receptor 2L13 protein (158-194AA)

Storage

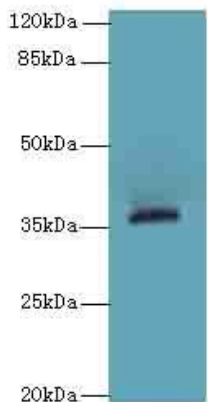
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Olfactory receptor 2L13, Olfactory receptor 2L14, OR2L13, OR2L14

Application

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



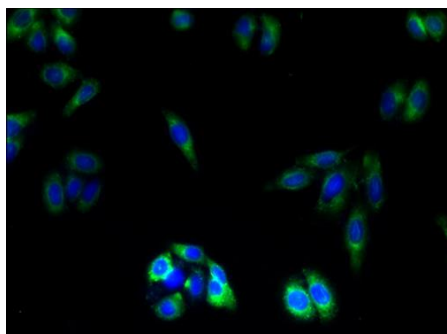
Western blot

All lanes: OR2L13 Polyclonal Antibody at 6ug/ml+Mouse liver tissue

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 36 kDa

Observed band size: 36 kDa



Immunofluorescent analysis of HepG2 cells using OR2L13 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)