
OR2L13 Polyclonal Antibody, FITC Conjugated

(Catalog #A66352)

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, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

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; Recommended dilution: ELISA 1:100-1:500
