

## TAS2R14 Polyclonal Antibody

(Catalog A67414)

### Background

Receptor that may play a role in the perception of bitterness and is gustducin-linked. May play a role in sensing the chemical composition of the gastrointestinal content. The activity of this receptor may stimulate alpha gustducin, mediate PLC-beta-2 activation and lead to the gating of TRPM5.

### Description

TAS2R14 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9NYV8

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Taste receptor type 2 member 14 protein (283-317AA)

### Storage

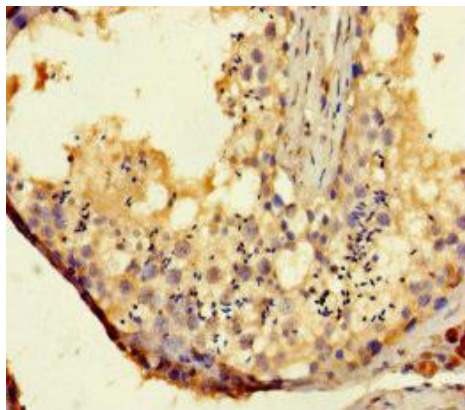
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Taste receptor type 2 member 14, T2R14, Taste receptor family B member 1, TRB1, TAS2R14

### Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human testis tissue using TAS2R14 Polyclonal Antibody at dilution of 1:100