

OPN5 Polyclonal Antibody

(Catalog # A65386)

Background

OPN5 (Opsin 5) is a member of the guanine nucleotide-binding protein (G protein)-coupled receptor superfamily. Expressed in the eye, brain, testes, and spinal cord, this gene belongs to the seven-exon subfamily of mammalian opsin genes that includes peropsin (RRH) and retinal G protein-coupled receptor (RGR). Similar to other seven-exon opsin genes, this family member may encode a protein with photoisomerase activity. Alternative splicing results in multiple transcript variants.

Description

OPN5 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

Q6U736

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Opsin-5 protein (310-353AA)

Storage

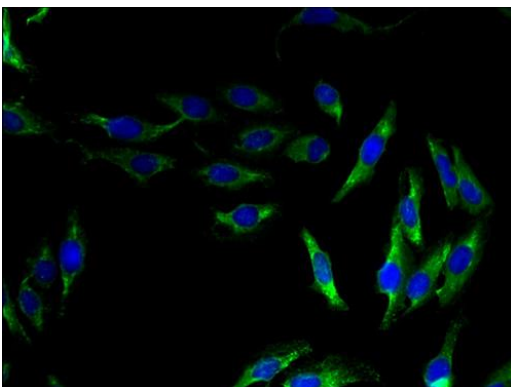
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

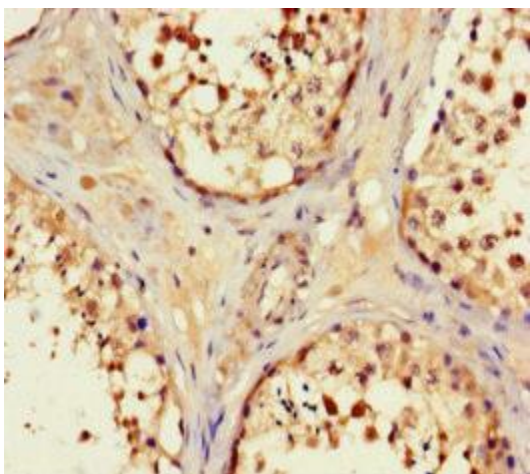
Opsin-5, G-protein coupled receptor 136, G-protein coupled receptor PGR12, Neuropsin, Transmembrane protein 13, OPN5, GPR136, PGR12, TMEM13

Application

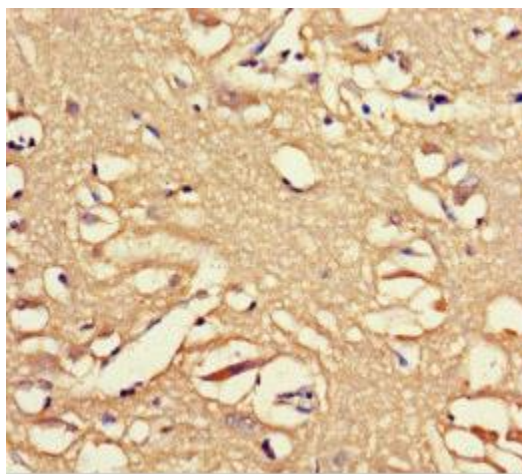
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



All lanes: FGF21 Polyclonal Antibody at 2ug/ml+A549 whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/15000 dilution
Predicted band size:
23 kDa Observed band size: 23 k



Immunohistochemistry of paraffin-embedded human testis tissue using OPN5 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human brain tissue using OPN5 Polyclonal Antibody at dilution of 1:100