
MS4A8 Polyclonal Antibody

(Catalog #A67798)

Background

May be involved in signal transduction as a component of a multimeric receptor complex.

Description

MS4A8 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

Q9BY19

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Membrane-spanning 4-domains subfamily A member 8 protein (20-74AA)

Storage

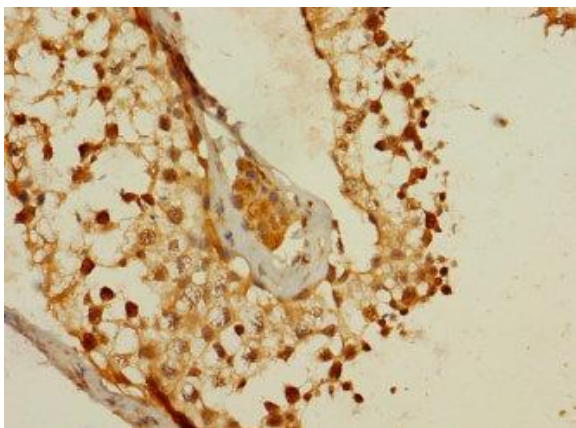
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

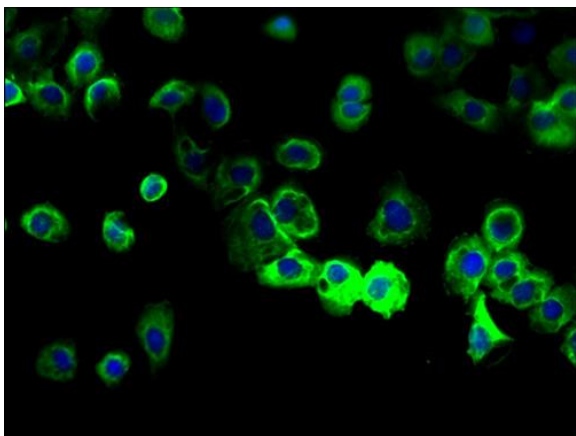
Membrane-spanning 4-domains subfamily A member 8, Four-span transmembrane protein 4, Membrane-spanning 4-domains subfamily A member 8B, MS4A8, 4SPAN4, MS4A8B

Application

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



Immunohistochemistry of paraffin-embedded human testis tissue using MS4A8 Polyclonal Antibody at dilution 1:100



Immunofluorescent analysis of A431 cells using MS4A8 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)