

# GGPS1 Polyclonal Antibody

(Catalog # A50241)

## Background

Catalyzes the trans-addition of the three molecules of IPP onto DMAPP to form geranylgeranyl pyrophosphate, an important precursor of carotenoids and geranylated proteins.

## Description

GGPS1 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

O95749

## Purification

>95%, Protein G purified

## Immunogen

Recombinant Human Geranylgeranyl pyrophosphate synthase protein (1-300AA)

## Storage

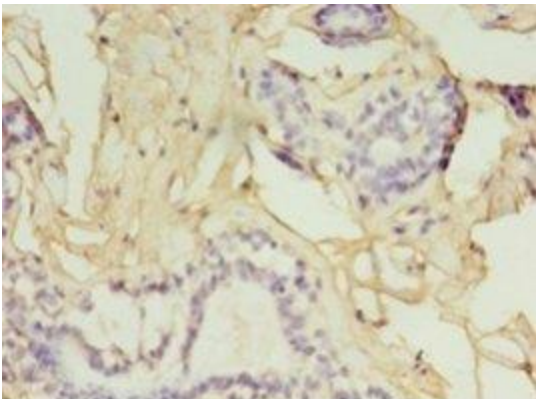
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

## Alternative Names

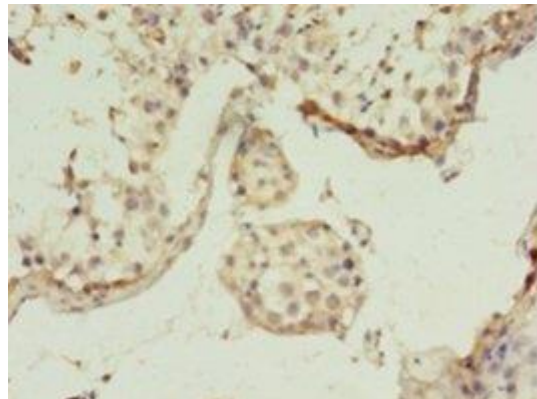
GGPP synthase, GGPPSase, (2E, 6E)-farnesyl diphosphate synthase, Dimethylallyltranstransferase, Farnesyl diphosphate synthase, Farnesyltranstransferase, Geranylgeranyl diphosphate synthase, Geranyltranstransferase, GGPS1

## Application

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



Immunohistochemistry of paraffin-embedded human breast cancer at dilution 1:100



Immunohistochemistry of paraffin-embedded human testis tissue at dilution 1:100