

PIGZ Polyclonal Antibody

(Catalog A65878)

Background

Mannosyltransferase involved in glycosylphosphatidylinositol-anchor biosynthesis. Transfers a fourth mannose to some trimannosyl-GPIs during GPI precursor assembly. The presence of a fourth mannose in GPI is facultative and only scarcely detected, suggesting that it only exists in some tissues.

Description

PIGZ 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

Q86VD9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human GPI mannosyltransferase 4 protein (1-130AA)

Storage

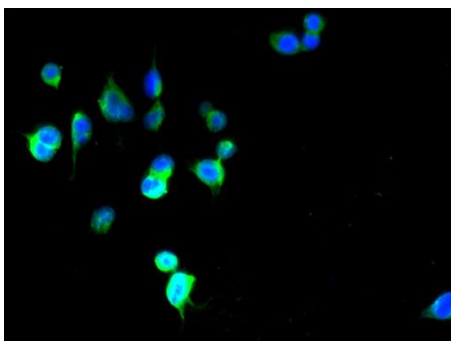
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

GPI mannosyltransferase 4, GPI mannosyltransferase IV, GPI-MT-IV, Phosphatidylinositol-glycan biosynthesis class Z protein, PIG-Z, SMP3 homolog, hSMP3, PIGZ, SMP3

Application

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



Immunofluorescent analysis of 293T cells using PIGZ Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)