
Gp2 Polyclonal Antibody

(Catalog # A67678)

Background

apical plasma membrane, extracellular exosome, antigen binding, antigen transcytosis by M cells in mucosal-associated lymphoid tissue

Description

Gp2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

Q9D733

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse Pancreatic secretory granule membrane major glycoprotein GP2 protein (22-531AA)

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

Pancreatic secretory granule membrane major glycoprotein GP2, Pancreatic zymogen granule membrane protein GP-2, Gp2

Application

ELISA
