

Echelon Biosciences Inc.

Mouse Anti-PIP_n Antibody

Z-P999

Support: echelon@echelon-inc.com

Description:

Mouse monoclonal antibody targeting PIPs

Applications:

ELISA - 0.5 ug/mL

Lipid-Protein Overlay - 1.0 ug/mL

Other in vitro and cellular applications are possible using this antibody, but have not been verified by Echelon Biosciences.

Properties:

Form - liquid

Storage instructions - Store at 4 °C for up to 30 days. Aliquot and store at -20 or -70 °C if longer storage is necessary. Avoid repeated freeze/thaw cycles.

Storage buffer - PBS, pH 7.4

Concentration - 1.0 mg/mL

Purity - affinity purified

Immunogen - synthetic PIP

Clonality - monoclonal

Isotype - IgG2b

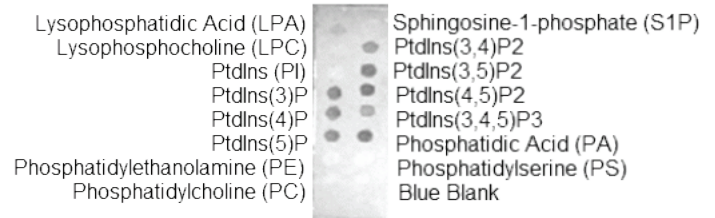
Specificity:

Z-P999 reacts with all seven phosphoinositides, and in some applications, phosphatidylinositol.

Background:

Phosphoinositides (PIPns) are minor components of cellular membranes but are integral signaling molecules for cellular communication.

Data: PIP Strip (Echelon P-6001)



Z-P999 was run on a PIP Strip at 1 ug/mL in PBS-T (0.1% v/v Tween-20) buffer. The membrane was not pre-blocked as the antibody does not work well under these conditions. The secondary antibody was incubated using PBS-T and developed using chemiluminescence.

Related Products:

Products	Catalog Number
Assays, Lipids, and Enzymes	
PIP3 Mass ELISA	K-2500s
PI3-Kinase Activity ELISA	K-1000s
PI3Ka enzyme, active	E-2000
Lipids and Antibodies	
PI(3,4,5)P3	P-3908, P-3916
PI(4,5)P2	P-4508, P-4516
PI(4)P	P-4008, P-4016
PI(3)P	P-3008, P-3016
Anti-PI(3,4,5)P3	Z-P345 (purified IgM)
Anti-PI(4,5)P2	Z-P045 (purified IgM)

Technical Data Sheet Rev. 2, 07-30-18 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.com