

Echelon Biosciences Inc.

Anti-Phosphatidylethanol Antibody

Z-PETH

Support: echelon@echelon-inc.com

Description:

Recombinant bivalent human monoclonal F(ab)₂ targeting phosphatidylethanol (PEth)

Z-PETH is a F(ab) dimer containing a heavy chain C-terminal dHLX dimerization domain (HuCAL®).

Applications:

ELISA - 0.5 ug/mL

IF - 2.5 ug/mL¹

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 – 3x PBS, pH 7.4

Concentration – 0.25-1.0 mg/mL. See product label.

Purity – affinity purified

Immunogen – synthetic PEth

Clonality – monoclonal, clone 367

Isotype – IgG

Molecular Weight – 115 kDa

Tag: Myc, Strep or FLAG, His6. This is lot specific. Please inquire.

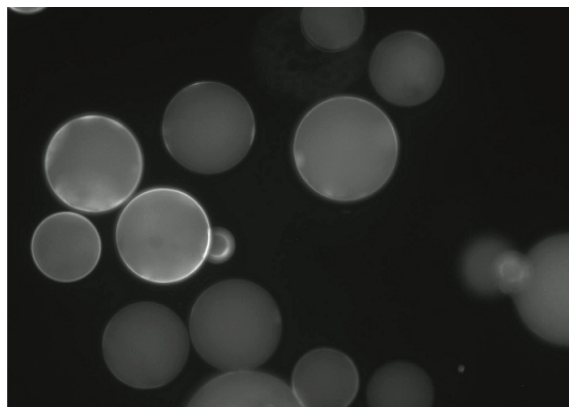
Specificity:

Z-PETH reacts primarily with phosphatidylethanol (of synthetic or natural origin), and demonstrates low cross-reactivity with other lipids and phospholipids including phosphatidylcholine, phosphatidylethanolamine, phosphatidic acid, phosphatidylserine, phosphatidylglycerol, sphingomyelin, and cholesterol.

Background:

PEth is formed in the presence of ethanol via the enzymatic action of phospholipase D on phosphatidylcholine and other phospholipids. Circulating levels of PEth can be used as a biomarker to measure alcohol consumption.

Data: Immunofluorescence



FITC labeled Z-PETH bound to PEth coated agarose beads (Echelon Cat # P-BPEth).

References:

1. Esteban-Pretel, G., M. P. Marin, et al. (2012). "Polyphosphoinositide Metabolism and Golgi Complex Morphology in Hippocampal Neurons in Primary Culture is Altered by Chronic Ethanol Exposure." Alcohol and Alcoholism.

Related Products:

Products	Catalog Number
Assays, Reagents, and Lipids	
PEth ELISA	K-5000
PEth	L-6017
Biotinylated PEth	L-60B16
BODIBY-FL PEth	L-60F18
PEth Beads	P-BPEth

Technical Data Sheet Rev. 5, 06-06-23 - 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