

Mouse Anti-PI(3,4)P₂ Antibody

Z-P034

Support: echelon@echelon-inc.com

Description:

Mouse monoclonal antibody targeting PI(3,4)P₂

Applications:

IF – 5 µg/mL
Lipid-Protein Overlay – 0.5 µg/mL
ELISA – 0.5 µg/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 °C if longer storage is necessary. Avoid repeated freeze/thaw cycles.

Storage buffer – PBS, pH 7.4

Concentration – see product label

Purity – affinity purified

Immunogen – synthetic PI(3,4)P₂

Clonality – monoclonal

Specificity:

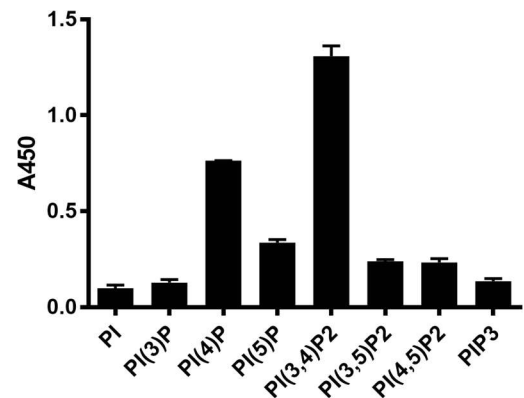
Z-P034 reacts primarily with the head group of the indicated phosphoinositide (of synthetic or natural origin) and demonstrates low cross-reactivity with other phosphoinositide or phospholipid depending on the assay format.

Background:

Phosphatidylinositol 3,4-bisphosphate (PI(3,4)P₂) is produced in stimulated cells by the action of phosphatidylinositol 3-kinases (PI3Ks) and/or lipid phosphatases (SHIP,

TPTE, etc.). PI(3,4)P₂ binds the pleckstrin homology (PH)-domain of AKT/PKB leading to cell survival and it also has an important role in podosome formation near focal adhesions.

Data: ELISA (specificity using Echelon H-6300)



References:

1. Rajala RV, Rajala A, Morris AJ, Anderson RE (2014) Phosphoinositides: minor lipids make a major impact on photoreceptor cell functions. *Sci Rep.* ;4:5463.
2. Chiang HC, Wang L, Xie Z, Yau A, Zhong Y. (2010) PI3 kinase signaling is involved in Abeta-induced memory loss in Drosophila. *Proc Natl Acad Sci U S A.* ;107(15):7060-5. PubMed PMID: 20351282.

Related Products:

Products	Catalog Number
Assays and Reagents	
PI(3,4)P ₂ PolyPIPosomes	Y-P034
PI(3,4)P ₂ Beads	P-B034a
PI(3,4)P ₂ Mass ELISA	K-3800
Lipids and Antibodies	
PI(3,4)P	P-3408, P-3416
Anti-PI(3,4)P ₂	Z-P034b