

Echelon Biosciences Inc.

Mouse Biotinylated Anti-PI(3,4)P₂ Antibody

Z-B034

Support: echelon@echelon-inc.com

Description:

Mouse monoclonal biotinylated antibody against PI(3,4)P₂

Applications:

ELISA - 0.5 ug/ml
Lipid-protein overlay - 0.5 ug/ml
AlphaScreen - 0.25-1.0 ug/ml
FP - 250 ug/ml
IF/ICC - 5.0 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.
Concentration - 0.5 mg/mL
Purity - affinity purified
Immunogen - synthetic PI(3,4)P₂
Clonality - monoclonal
Isotype - IgG3

Specificity:

Z-B034 reacts primarily with the head group of the indicated phosphoinositide, and demonstrates low cross-reactivity with other phosphoinositides or phospholipids 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.

References:

1. Chiang, H. C., L. Wang, et al. (2010). PI3 kinase signaling is involved in Abeta-induced memory loss in Drosophila. PNAS 107(15): 7060-5.
2. Bae, Y. H., Z. Ding, et al. (2010). Profilin1 regulates PI(3,4)P₂ and lamellipodin accumulation at the leading edge thus influencing motility of MDA-MB-231 cells. Proceedings of the National Academy of Sciences.

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-P034 Z-P034b

Technical Data Sheet Rev. 3, 06-20-18 - 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