

Echelon Biosciences Inc.

Mouse HRP-conjugated Anti-PI(3,4,5)P₃

Z-H345

Support: echelon@echelon-inc.com

Description:

Monoclonal mouse IgM antibody conjugated to HRP targeting PI(3,4,5)P₃

Applications:

IF - 5 ug/mL

ELISA - 20 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.

For long-term storage, aliquot and store at -20 °C

Avoid repeated freeze/thaw cycles.

Storage buffer - PBS + 0.01% thiomersal

Concentration - 2.7 mg/mL

Purity - affinity purified by anti-IgM chromatography

Immunogen - PI(3,4,5)P₃ conjugated to BSA

Clonality - monoclonal; RC6F8

Isotype - IgM

Specificity:

HRP Anti-PI(3,4,5)P₃ reacts primarily with the head group of the indicated phosphoinositide (of synthetic or natural origin), and demonstrates low cross-reactivity with other phosphoinositides or phospholipids depending on the assay format.

Background:

PI(3,4,5)P₃ is generated by phosphorylation of PI(4,5)P₂ by PI3 kinases. Production of PI(3,4,5)P₃ at the plasma membrane is known to enhance association of PH-domain containing proteins associated with multiple receptor signaling cascades.

References:

1. Silva A, Yunes JA, Cardoso BA, Martins LR, Jotta PY, Abecasis M, Nowill AE, Leslie NR, Cardoso AA, Barata JT. (2008) PTEN posttranslational inactivation and hyperactivation of the PI3K/Akt pathway sustain primary T cell leukemia viability. *J Clin Invest.* 118(11):3762-74.
2. Hama H, Torabinejad J, Prestwich GD, DeWald DB. (2004) Measurement and immunofluorescence of cellular phosphoinositides. *Methods Mol Biol.* 2;284:243-58.
3. Chen R, Kang VH, Chen J, Shope JC, Torabinejad J, DeWald DB, Prestwich GD. (2002) A monoclonal antibody to visualize PtdIns(3,4,5)P(3) in cells. *J Histochem Cytochem.* 50(5):697-708.

Related Products:

Products	Catalog Number
Assays, Lipids, and Enzymes	
PI3 Mass ELISA	K-2500s
PI3-Kinase Activity ELISA	K-1000s
SHIP2 Enzyme	E-1000
Lipids and Antibodies	
PI(3,4,5)P ₃	P-3908, P-3916
Anti-PI(3,4,5)P ₃	Z-P345b (purified IgG) Z-B345b (biotinylated IgG) Z-P345 (purified IgM)

Technical Data Sheet Rev. 5, 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