



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name

Anti-PI(4,5)P₂, in cell supernatant

Product Number:

Z-S045

Description:

Anti-phosphatidylinositol-4,5-bisphosphate [Anti-PtdIns(4,5)P₂]
IgM monoclonal antibody in cell supernatant.

Storage:

Store at 2-8°C for up to 30 days. Aliquot and freeze at -20°C if longer
storage is necessary. Do not freeze-thaw antibody.

Hybridoma Source:

2C11

Immunogen:

Synthetic PtdIns(4,5)P₂ (1) in liposomes.

Suggested Usage:

1:500 for dot blot. 1:250 for TLC overlay and ELISA. 1:50 for
cellular immunofluorescence and immunohistochemistry. Detailed
procedures are available from Echelon Biosciences Inc.

Antibody Specificity:

Shown to recognize PtdIns(4,5)P₂ in murine and human cells. Some
crossreactivity with PI4P, PI(3,4,5)P₃ and PI(3,5)P₂. Crossreactivity is
1% or less with other phosphatidylinositol phosphates or lipids.

Reference:

1. Thomas, C.L., Steel, J., Prestwich, G.D., and Schiavo, G.
Generation of Phosphatidylinositol-specific Antibodies and their
Characterization Biochemical Society Transactions (1999) 27 (4)
648-652

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

TDS Z-S045 Rev: 3 (07/01/03)

**Use this
Product with:**

D-myo-phosphatidylinositol 4,5-bisphosphates (PtdIns(4,5)P₂)
including:

- D(+)-sn-1,1-di-*O*-butanoylglyceryl,3-*O*-phospho linked (Product P-4504)
- D(+)-sn-1,1-di-*O*-octanoylglyceryl,3-*O*-phospho linked (Product P-4508)
- D(+)-sn-1,1-di-*O*-hexadecanoylglyceryl,3-*O*-phospho linked (Product P-4516)
- NBD6 PtdIns(4,5)P₂ Fluorescent Probe

D(+)-sn-1-*O*-[6'-[[6-[(7-nitrobenz-2-oxa-1,3-diazol-4-yl)amino]hexanoyl]amino]-hexanoyl]-2-*O*-hexanoylglyceryl *D-myo*-phosphatidylinositol 4,5-bisphosphate (Product C45N6)

**Related Products
In Development:**

Anti-phosphatidylinositol-4,5-bisphosphate [Anti-PtdIns(4,5)P₂]
purified from ascites (Product No. Z-P45)

Anti-phosphatidylinositol-4,5-bisphosphate [Anti-PtdIns(4,5)P₂] in
ascites (Product No. Z-S45)

Anti-phosphatidylinositol-4,5-bisphosphate [Anti-PtdIns(4,5)P₂]
enzyme or fluorescent conjugates

A full line of antibodies specific to phosphatidylinositol phosphates