



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

Certificate of Analysis

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Custom Lipid Array Strip

Catalog number: P-6111-P-3016

Number of Strips: 30

Lipid Spotted: Phosphatidylinositol 3-phosphate diC16, catalog # P-3016

Quality Control Protein (if applicable*): PI(3)P Grip (P40PX), Catalog # G-0302

Storage: Store at 2-8 °C. Product is moisture and light sensitive.

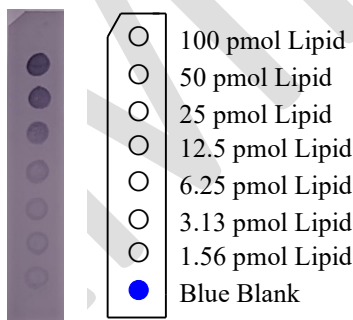
Product Description:

1.25 x 6 cm hydrophobic membrane that has been spotted with 7 concentrations of Lipid and a no lipid control (100 pmol, 50, 25, 12.5, 6.25, 3.13, 1.56, 0). These membranes can be used to determine lipid-protein interactions, through a simple protein-lipid overlay experiment.

Format:

The membrane has a diagonal cut on its top left corner and is spotted with Xylene Cyanol FF (blue) at the bottom to assist in orientation of the strip. Ponceau S staining (pink)** was added to the lipid spots. See template below for location of lipids.

Membrane Template and Quality Data:



Suggested Usage:

Visit our website www.echelon-inc.com. At the bottom of each product's webpage is our general Protocol "Protocol_Strip_Array" for use with product numbers: P-6001, P-6100, P-6002, P-6003, S-6000, and S-6001. Also please refer to our FAQ "Frequently Asked Questions" document.

References:

1. Dowler S, Currie RA, Downes PC, Alessi DR. DAPP1: a dual adaptor for phosphotyrosine and 3-phosphoinositides. *Biochemical Society J.* 342, 7-12 (1999)
2. Dowler, S., Kular, G., and Alessi, R.D., Protein lipid overlay assay, *Sci STKE*, 2002. April 23; 2002 (129). p16.
3. Ferguson CG, James RD, Bigman CS, Shepard DA, Abdiche Y, Katsamba PS, Myszka DG, Prestwich GD. Phosphoinositide-containing polymerized liposomes: stable membrane-mimetic vesicles for protein-lipid binding analysis. *Bioconjug Chem.* 2005 Nov – Dec; 16(6): 1475-83.
4. Busse RA, ScaciocA, Hernandez JM, Krick R, Stephan M, Janshoff A, Thumm M, Kuhnel K. Qualitative and quantitative characterization of protein-phosphoinositide interactions with liposome-based methods. *Autophagy*. 2013 May 1;9(5):770-7.

*Quality Control with a lipid binding protein will not be run for lipids without an internal binding protein.

**Final concentration of 0.002% (w/v) Ponceau S was added for accuracy during spotting.

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.