

# Echelon Biosciences Inc.

## Membrane Lipid Arrays™

P-6003

### Description

Membrane Lipid Arrays™ are 4 cm x 5 cm hydrophobic membranes that have been spotted with a concentration gradient of eight different lipids. These membranes can be used to determine lipid-protein interactions, through a simple protein-lipid overlay experiment. This allows researchers a convenient way to determine if their protein of interest interacts with one or more of the bound lipids.

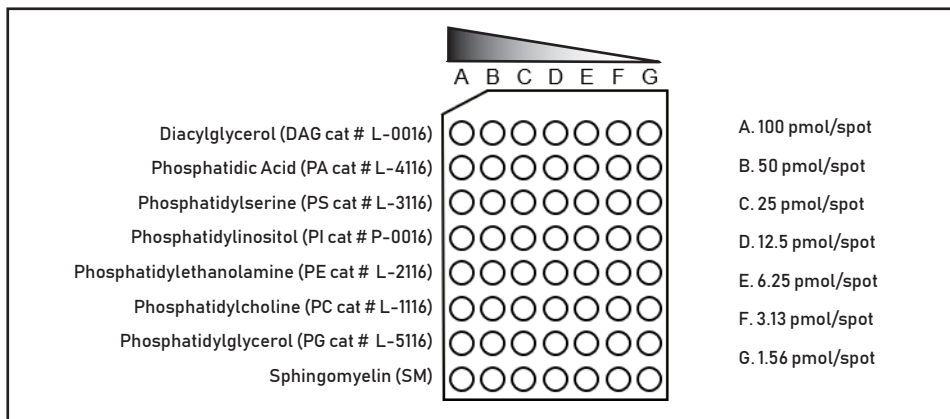
### Storage

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

### Format

The membrane has a diagonal cut on its top left corner and Ponceau S staining (pink) was added to the lipid spots to assist in orientation of the array. See template below for location of lipids. All of the lipids are long chain (diC16) highly pure natural or synthetic analogs. For more information, on the lipids spotted on the membrane, please visit our website and search the catalog numbers provided in the figure below.

### Membrane Template



### Suggested Use

Visit our website [www.echelon-inc.com](http://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, P-6111, 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. Dibble CF, Horst JA, Malone MH, Park K, Temple B, Cheeseman H, Barbaro JR, Johnson GL, Bencharit S. Defining the functional domain or programmed cell death 10 through its interactions with phosphatidylinositol-3,4,5-triphosphate. *PLoS One.* 2012 Jul 23;5(7):e11740.

### Warranty and Disclaimer:

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. This product has a quality warranty of 6 months from the date of receipt when stored properly. We welcome inquiries about licensing the use of our trademarks and technologies at [busdev@echelon-inc.com](mailto:busdev@echelon-inc.com).

Technical Data Sheet Rev. 3, 08-03-18 - 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](http://Echelon-inc.com)