

Echelon Biosciences Inc.

Mega Lipid Strips™

P-6005

Description

Mega Lipid Strips™ are 2.5 cm x 6 cm hydrophobic membranes that have been spotted with 23 biologically relevant lipids at 100 pmol per spot. 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

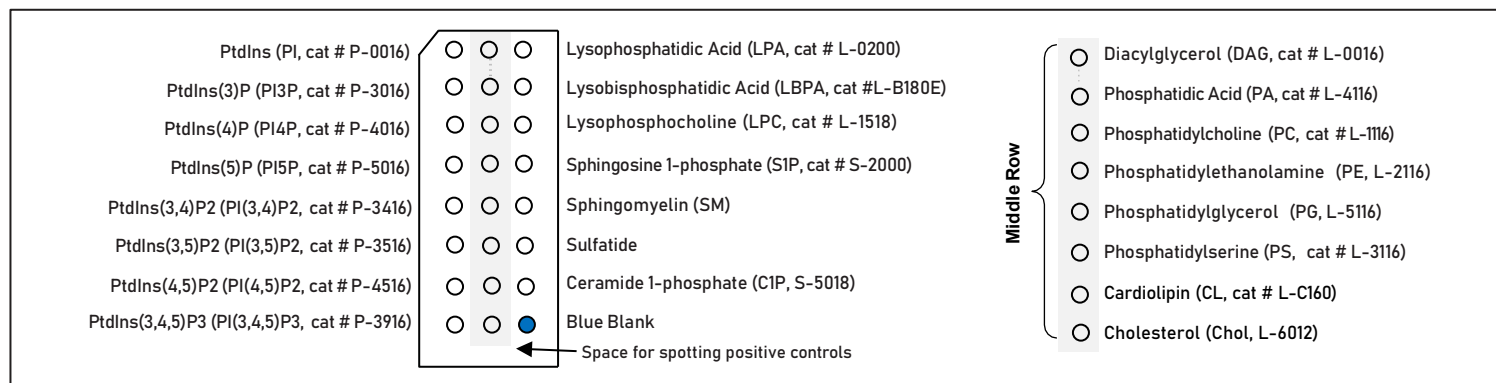
Store at 2-8 °C. Product is moisture and light sensitive.

Format

The membrane has a diagonal cut on its top left corner and is spotted with Xylene Cyanol FF* (blue) in the bottom right blank spot to assist in orientation of the strip. Ponceau S staining (pink) was added to the lipid spots. Some space is provided at the bottom of each Mega Lipid Strip for additional control proteins or antibodies to be spotted by the customer. See template below for location of lipids. All the lipids are long chain (\geq C16) 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.

*Xylene Cyanol FF shows fluorescence around 615 nm and may be visible when using fluorescence-based imaging such as Licor.

Membrane Template



Visit website www.echelon-inc.com for protocols and frequently asked questions.

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

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.

Technical Data Sheet Rev. 1, 03-08-24 - 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