

Echelon Biosciences Inc.

Sphingo Strips™

S-6000

Description

Sphingo Strips™ are 2 cm x 6 cm hydrophobic membranes that have been spotted with 15 different 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 PIP Strip for additional control proteins or antibodies to be spotted by the customer. See template below for location of lipids. All of 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

Sphingosine (cat # S-1000)	<input type="radio"/>	<input type="radio"/>	Monosialoganglioside-GM1
Sphingosine 1-phosphate (S1P, cat # S-2000)	<input type="radio"/>	<input type="radio"/>	Disialoganglioside-GD3
Phytosphingosine	<input type="radio"/>	<input type="radio"/>	Sulfatide
Ceramide	<input type="radio"/>	<input type="radio"/>	Psychosine
Sphingomyelin (SM)	<input type="radio"/>	<input type="radio"/>	Cholesterol (cat # L-6012)
Sphingosylphosphorylcholine (SPC)	<input type="radio"/>	<input type="radio"/>	Lysophosphocholine (LPC, cat # L-1518)
Lysophosphatidic Acid (LPA, cat # L-0200)	<input type="radio"/>	<input type="radio"/>	Phosphatidylcholine (PC, cat # L-1116)
Myriocin	<input type="radio"/>	<input checked="" type="radio"/>	Blue Blank

Suggested Use

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, P-6111, S-6000, and S-6001. Also please refer to our FAQ "Frequently Asked Questions" document.

References

1. Dowler, S., Currie, R.A., Downes, P.C., and Alessi, D.R. 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, L6. (2002).
3. Rodriguez-Asiain A, Ruiz-Babot G, Romero W, Cubi R, Erazo T, Biondi RM, Bavascas JR, Aquilera J, Gomez N, Gil C, Claro E, Lizcano JM. Brain specific kinase-1 BRSK1/SAD-B associates with lipid rafts: modulation of kinase activity by lipid environment. *Biochem Biophys.* 2011 Dec.
4. Sawa CG, Fernandes da Costa SP, Bokori-Brown M, Naylor CE, Cole AR, Moss DS, Titball RW, Basak AK. Molecular architecture and functional analysis of NetB, a pore-forming toxin from *Clostridium perfringens*. *J Biol Chem.* 2013 Feb.

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. 9, 08-03-18 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Ph: 866-588-0455
Fax: 801-588-0497

Echelon-inc.com