



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:** SpingoStrips™

**Product Number:** S-6000

**General Description:** SpingoStrips™ are 2 x 6 cm hydrophobic membranes that have been spotted with 15 different biologically active 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 Ponceau S\* staining to assist in orientation of the strip. See template below for location of lipids.

**Membrane Template:**

Sphingosine	○	○	Monosialoganglioside-GM1
Sphingosine-1-phosphate (S1P)	○	○	Disialoganglioside-GD3
Phytosphingosine	○	○	3-sulfogalactosylceramide (Sulfatide)
Ceramide	○	○	Psychosine
Sphingomyelin	○	○	Cholesterol
Sphingosylphosphorylcholine (SPC)	○	○	Lysophosphocholine (LPC)
Lysophosphatidic Acid (LPA)	○	○	Phosphatidylcholine (PC)
Myriosine	○	○	Blank

**Suggested Usage:** See support protocol “basic Protocol for PIP strips™, PIP Array™, PIP Microstrips™, and SpingoStrips™,” on our website [www.echelon-inc.com](http://www.echelon-inc.com)

**References:** 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).  
Deak, M., Casamayor, A., Currie, R.A., Downes, P., Alessi, Dowler, S., Kular, G., and Alessi, R.D., Protein lipid overlay assay, *Sci STKE*, 2002, L6. (2002).

\*Final concentration of 0.1% (v/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](mailto:busdev@echelon-inc.com).