



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:** PIP Array™

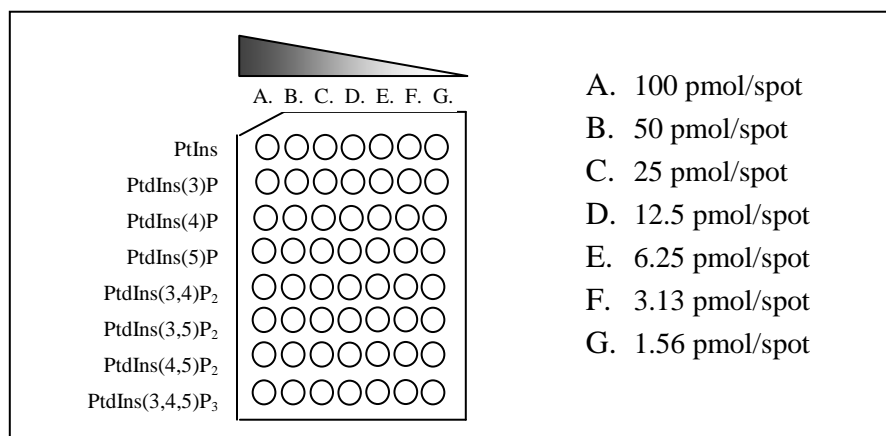
**Product Number:** P-6100

**General Description:** PIP Array™ are 4 x 5 cm hydrophobic membranes that have been spotted with a concentration gradient of all eight phosphoinositides. 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 and the degree of binding.

**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:



**Suggested Usage:** See support protocol “basic Protocol for PIP strips™, PIP Array™, PIP Microstrips™, and SphingoStrips™” 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).  
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.