



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.

## Cova PIP™ Screening Plate

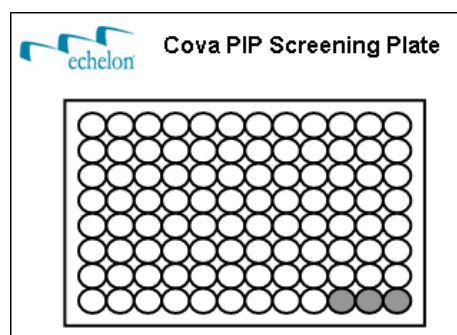
*Phosphoinositide-coated microtiter plates*

**Product Number: H-6200**

**Description:** Transparent 96-well (12 x 8 strip-well) polystyrene microtiter plate containing 20 pmols of an individual diC<sub>6</sub>-modified PIP<sub>n</sub> per well. Three wells are left without lipid, but still blocked, for a negative control. The coated volume is 100 µL; and each plate is pre-blocked and ready for addition of protein. See figure.

Lipids Available:

Cat #	Lipid
H-6201	PI
H-6203	PI(3)P
H-6204	PI(4)P
H-6205	PI(5)P
H-6234	PI(3,4)P <sub>2</sub>
H-6235	PI(3,5)P <sub>2</sub>
H-6245	PI(4,5)P <sub>2</sub>
H-6239	PI(3,4,5)P <sub>3</sub>



**Storage:** Store at 2-8°C for up to 1 year after date of receipt. Product is sensitive to moisture.

**Suggested Use:** Determination of protein interaction with PIP<sub>n</sub>s via solid phase binding assays.

**QA/Product Testing:** Cova PIP™ Screening Plates are tested for deposition of phosphoinositides and the expected binding-selectivity of certain PH-domain proteins (PIP Grip™ product line) or purified anti-phosphoinositide antibodies (Z-P### product line). For additional information, visit our web site ([www.echelon-inc.com](http://www.echelon-inc.com)) or contact our technical service department at 1-866-588-0455.

### Related Products:

- Cova PIP Specificity Plates (H-6300)
- Cova PIP Sensitivity Plates (H-6400)
- Cova PIP Array Plates (H-6500)
- Lipid Coated Beads (P-Bxxx)
- Polymerized liposomes, PolyPIPosomes (Y-Pxxx)
- PIP Strips (P-6001), PIP-Arrays (P-6100).