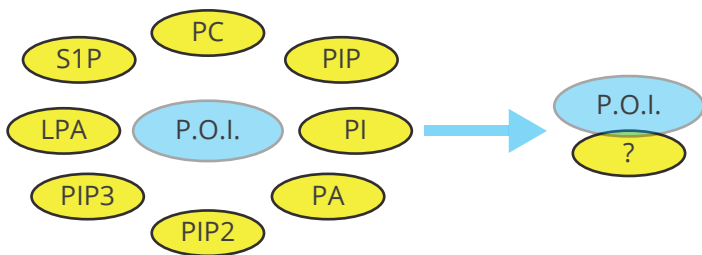


## Does your protein of interest bind lipids?



Lipids are fundamental to cell biology, both as the building blocks for biological membranes and as intracellular signaling molecules. A less appreciated, but equally important feature of lipids is their ability to bind and influence the structure and function of proteins. While these interactions are critical, studying them can be challenging.

**Echelon Biosciences** has developed a suite of lipid-protein interaction tools that lower the barriers for interrogating lipid-protein binding partners.

## Lipid Strips

**PIP Strips** (Cat # P-6001) are 2 x 6 cm hydrophobic membranes that have been spotted with 100 pmol of all eight phosphoinositides and seven other biologically important lipids. These strips can be used to determine if a protein of interest interacts with one or more of these lipids in a simple protein-lipid overlay assay.

Also available:

**PIP Array** (P-6100), **Membrane Lipid Strips** (P-6002), **Membrane Lipid Array** (P-6003), **Sphingo Strips** (S-6000), **Sphingo Array** (S-6001), **Custom PIP Array** - contact us

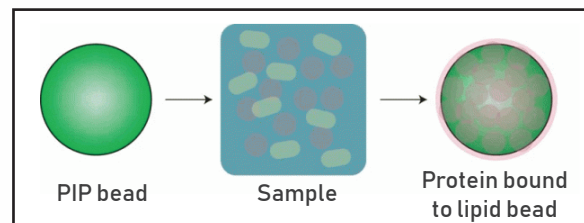
## PIP Strips™ (Cat # P-6001)

Lysophosphatidic Acid	<input type="radio"/>	<input type="radio"/>	Sphingosine-1-phosphate
Lysophosphocholine	<input type="radio"/>	<input type="radio"/>	PtdIns(3,4)P2
PtdIns	<input type="radio"/>	<input type="radio"/>	PtdIns(3,5)P2
PtdIns(3)P	<input type="radio"/>	<input type="radio"/>	PtdIns(4,5)P2
PtdIns(4)P	<input type="radio"/>	<input type="radio"/>	PtdIns(3,4,5)P3
PtdIns(5)P	<input type="radio"/>	<input type="radio"/>	Phosphatidic Acid
Phosphatidylethanolamine	<input type="radio"/>	<input type="radio"/>	Phosphatidylserine
Phosphatidylcholine	<input type="radio"/>	<input checked="" type="radio"/>	Blue Blank

## Lipid Beads

**Lipid Beads** are ideal for use in protein pull-down experiments to identify and characterize lipid binding proteins. The beads allow lipid binding proteins to be isolated from cell lysates or mixtures of in vitro translated peptides, which can then be analysed by SDS-PAGE, Western blot, or mass spectrometry. See the table for available Lipid Beads.

Visit [echelon-inc.com](http://echelon-inc.com) for more Lipid Beads. Contact us for **custom** bead conjugations.



Lipid Coated Beads from Echelon			
Lipid	Cat #	Lipid	Cat #
Control	P-B000	PI(3,5)P2	P-B035a
PtdIns	P-B001	PI(4,5)P2	P-B045a
PI(3)P	P-B003a	PI(3,4,5)P3	P-B345a
PI(4)P	P-B004a	PA	P-B0PA
PI(5)P	P-B005a	PC	P-B0PC
PI(3,4)P2	P-B034a	PS	P-B0PS

