

Lipid Coated Beads

Support: echelon@echelon-inc.com

Description: Lipid Beads™ are designed for protein pull-down experiments. Possible uses are to test for lipid binding in a solution of purified protein, in a cell lysate, or to test binding of radiolabeled *in vitro* translation products.

Storage: Store at 4°C. Do not freeze!

Lipid Bead	Catalog #
Control (No Lipid)	P-B000
PtdIns	P-B001
PI(3)P	P-B003a
PI(4)P	P-B004a
PI(5)P	P-B005a
PI(3,4)P2	P-B034a
PI(3,5)P2	P-B035a
PI(4,5)P2	P-B045a
PI(3,4,5)P3	P-B345a
Cardiolipin	P-BCLP
Monolysocardiolipin	P-BMCL
Ceramide 1-phosphate	P-BC1P
Ceramide	P-BCer
Glucosylceramide	P-B0GC
LBPA	P-BLBPA
Lactosyl Ceramide	P-B0LC

Lipid Bead	Catalog #
Phosphatidic Acid	P-B0PA
Phosphatidylcholine	P-B0PC
Phosphatidylethanolamine	P-B0PE
Phosphatidylglycerol	P-B0PG
Phosphatidylserine	P-B0PS
Sphingomyelin	P-B0SM
Phosphatidylethanol	P-BPEth
Sulfatide	P-Bsulf
Cholesterol	P-BChl
Lysophosphatidic Acid	L-6101
Ins(1,3,4,5)P4	Q-B1345
Sphingosine	S-6100
Sphingosine 1-phosphate	S-6110
PIP Beads Sample Pack	P-B00S
Sphingo Beads Sample Pack	P-B00SS

Lipid Bead™ Assay Notes

1. Lipid Coated Beads are available in 1 mL and 0.1 mL sizes and come as a 50% slurry in 1X PBS Buffer.
2. An order of 1 mL comes with 200 µL of Control Beads. Control Beads are also available for purchase and their use is suggested for analysis.
3. Store at 2-8°C. Product is temperature and light sensitive. Do not freeze.
4. Centrifuge the beads at 1,000 x g or lower. Do not centrifuge the beads at high speed as this can crush the beads.
5. There is a total of 10 nmoles of bound lipid per 1 mL of beads.
6. The beads range from 45 – 165 micrometers in diameter.
7. Use 50-100 µL of 50 % slurry lipid coated beads and control beads for each protein.
8. See table 1 for a list of positive control proteins.

Lipid Bead™ - Protein Pull-down Protocol

We provide the following protocol as a guide, and strongly encourage researchers to consult scientific literature and conduct optimization experiments to establish the most favorable conditions for their protein of interest.

1. Mix beads well (do not vortex). Transfer 50 – 100 µL of 50% bead slurry to a 0.5 – 1.5 mL tube.
2. Pellet beads by centrifugation at 1,000 x g or lower. Carefully remove supernatant to avoid losing beads.
3. Wash beads with 10X excess of the wash buffer. Incubate wash 1-2 min with gentle mixing. Repeat wash steps two times.
4. Carefully remove wash supernatant and resuspend the beads in binding buffer containing your protein.
 - a. For purified proteins, use 1-10 µg of protein diluted in enough binding buffer to form a 50% lipid bead slurry.
 - b. For cell lysates, extracts, subcellular fractions: use approximately 1 mg of total protein to 1 mL binding buffer. Add the entire 1 mL of prepared sample to your washed beads.
5. Incubate the protein-bead solution for 1-3 hours. The incubation can be done at room temperature or at 4° C, depending on the stability of your protein. Continuous motion is required to keep beads in suspension.
6. Pellet the beads and carefully remove supernatant. If needed, the supernatant can be used to monitor unbound protein.
7. Wash beads with 10X excess of the wash buffer. Incubate wash 1-2 min with gentle mixing. Repeat wash steps two to five times. If you choose to use BSA in your binding buffer, make sure it is washed out well before moving to the next step.
8. Following the last wash, elute bound proteins by adding an equal volume of 2X Laemmli sample buffer (or similar) to beads and heat to 70 °C for 10 minutes.
9. Following heating, pellet the beads and remove the supernatant. Store supernatant at 4°C until analysis. Discard beads.
10. Proteins can be separated by SDS-PAGE and analyzed by Coomassie blue staining of the gel, protein transfer and immunoblotting to detect proteins of interest, or autoradiography.

Suggested Buffer Systems

The buffers below are a good starting point for testing lipid protein interactions. Other buffers may show improved binding with your protein of interest. Protein stabilizers, such as 0.1-1% BSA, can stabilize some proteins and reduce non-specific lipid-protein interactions. Non-ionic detergents, such as Igepal and Tween 20, can reduce non-specific lipid interactions and reduce background binding to control beads.

Buffer System 1

Wash/Binding Buffer: 10 mM HEPES, pH 7.4, 150 mM NaCl, 0.25% Igepal

If the protein is sticking nonspecifically to Control Beads, try increasing the detergent concentration to 0.5% or higher.

Buffer System 2

Binding Buffer: 0.1-1% BSA in PBS with 0.05-0.1% Tween 20, pH 6.8-7.4

Wash Buffer: PBS with 0.05-1% tween 20, pH 6.8-7.4

Buffer System 3

Wash/Binding Buffer: PBS with 0.05-1% tween 20, pH 6.8-7.4

Table 1: Positive Control Proteins: Bead - protein pairs for use as positive controls in your experiments.

Lipid Bead	Bead Catalog #	Suggested Buffer System	Protein Catalog #	Lipid Binding Protein
PI(3)P Bead	P-B003a	2	G-0302	PI(3)P Grip
PI(4)P Bead	P-B004a	2	Z-P004	PI(4)P Antibody
PI(3,4)P2 Bead	P-B034a	1, 2	G-9901	MultiPIP Grip
PI(3,5)P2 Bead	P-B035a	2	G-3501	PI(3,5)P2 Grip
PI(4,5)P2 Bead	P-B045a	1, 2	G-4501	PI(4,5)P2 Grip
PI(3,4,5)P3 Bead	P-B345a	2	G-9901 G-3901	MultiPIP Grip PI(3,4,5)P3 Grip (Grp1-PH)
Phosphatidic Acid Bead	P-B0PA	2	G-9901	MultiPIP Grip
Monolysocardiolipin Bead	P-BMCL	3	G-CL01	Cardiolipin Grip
Cardiolipin Bead	P-BCLP	3	G-CL01	Cardiolipin Grip
Ins(1,3,4,5)P4 Bead	Q-B1345	1, 2	G-3901	PI(3,4,5)P3 Grip (Grp1-PH)
Lysophosphatidic Acid Bead	L-6101	2	Z-P200	Lysophosphatidic Acid Antibody

Table 2: Troubleshooting Guide

Issue	Suggestion
No binding	<ul style="list-style-type: none"> • Increase protein concentration • Increase incubation time • Add a protein stabilizer such as BSA • Protein has lost activity
More than one band	<ul style="list-style-type: none"> • Increase detergent concentration in wash/Binding Buffer • Increase number of washes • Add longer incubation steps per each wash
Binds Control bead	<ul style="list-style-type: none"> • Increase detergent concentration in wash/Binding Buffer • Increase number of washes • Add longer incubation steps per each wash

Preparation of Cell Lysate / Protein (Process may vary by cell type / source)

This is a general method for using cell lysate. Harsh lysis conditions should be avoided as they may interfere with lipid structure and binding. Non-ionic detergents such as Igepal, NP40, and Triton X-100 are less harsh than ionic detergents, non-denaturing, and used for solubilizing and purification of membrane protein complexes.

1. Isolate and centrifuge cells.
2. Lyse the pellet. (See below for lysis buffer suggestions)
3. Vortex and centrifuge to prepare clarified samples.

Lysis Buffer

Salts (10-20 mM HEPES, pH 7.4, 100-150 mM NaCl)

Non-ionic detergents* (such as 0.1 – 2% Igepal)

Divalent cations (0-10 mM EDTA and/or EGTA)

Protease inhibitors (complete mini protease tablet, Na orthovanadate, Na fluoride.)

Statement

The protein-pull down protocol has been tested with the beads and proteins contained in table 1. Protein-lipid interactions determined with Echelon's lipid coated bead products may vary from the binding interactions determined in other methods. We recommend researchers use a variety of methods to fully characterize the lipid binding preference of a particular protein. Echelon has several innovative products useful for determining protein-lipid interactions. These products include lipid overlay products such as PIP strips (P-6001), stabilized liposomes (PolyPIPosomes™ e.g. Y-P039) and PIP-Plates™ (e.g. H-6300).

References

1. Zhu Z, Chen J, Wang G, Elsherbini A, Zhong L, Jiang X, et al. Ceramide regulates interaction of Hsd17b4 with Pex5 and function of peroxisomes. *Biochimica et Biophysica Acta (BBA) - Molecular and Cell Biology of Lipids*. 2019.
2. Patel A, Kostyak J, Dangelmaier C, Badolia R, Bhavanasi D, Aslan JE, et al. ELMO1 deficiency enhances platelet function. *Blood Advances*. 2019;3(4):575-87.
3. Huang D, Cao L, Xiao L, Song J-x, Zhang Y-j, Zheng P, et al. Hypoxia induces actin cytoskeleton remodeling by regulating the binding of CAPZA1 to F-actin via PIP2 to drive EMT in hepatocellular carcinoma. *Cancer Letters*. 2019.
4. Hawse WF, Cattley RT. T cells transduce T-cell receptor signal strength by generating different phosphatidylinositols. *J Biol Chem*. 2019;294(13):4793-805. doi: 10.1074/jbc.RA118.006524. PubMed PMID: 30692200.
5. Fang C, Manes TD, Liu L, Liu K, Qin L, Li G, et al. ZFYVE21 is a complement-induced Rab5 effector that activates non-canonical NF-κB via phosphoinositide remodeling of endosomes. *Nature Communications*. 2019;10(1):2247.
6. Dudley LJ, Cabodevilla AG, Makar AN, Sztacho M, Michelberger T, Marsh JA, et al. Intrinsic lipid binding activity of ATG16L1 supports efficient membrane anchoring and autophagy. *The EMBO Journal*. 2019:e100554. doi: 10.15252/em-bj.2018100554.
7. Ventimiglia LN, Cuesta-Geijo MA, Martinelli N, Caballe A, Macheboeuf P, Miguet N, et al. CC2D1B Coordinates ESCRT-III Activity during the Mitotic Reformation of the Nuclear Envelope. *Developmental Cell*. 2018;47(5):547-63.e6.
8. Ulicna L, Kalendova A, Kalasova I, Vacik T, Hozák P. PIP2 epigenetically represses rRNA genes transcription interacting with PHF8. *Biochimica et Biophysica Acta (BBA) - Molecular and Cell Biology of Lipids*. 2018;1863(3):266-75.
9. Stein CS, Jadia P, Zhang X, McLendon JM, Abouassaly GM, Witmer NH, et al. Mitoregulin: A lncRNA-Encoded Micro-protein that Supports Mitochondrial Supercomplexes and Respiratory Efficiency. *Cell Reports*. 2018;23(13):3710-20.e8.
10. Nair S, Sng J, Boddupalli CS, Seckinger A, Chesi M, Fulciniti M, et al. Antigen-mediated regulation in monoclonal gammopathies and myeloma. *JCI insight*. 2018;3(8).