

## Protocol for immunocytochemistry (ICC) with lipid antibodies

This protocol is recommended as a general guide for immunostaining of cells with antibodies for detecting lipids of interest. While this procedure may be suitable for staining proteins in some systems, we do not recommend it as a starting point. Please see our alternative protocols for ICC with protein antibodies.

**Immunostaining of cellular lipids requires more specific procedures than are provided in general ICC protocols owing to the different cellular compartments in which they are found and how those compartments are affected by fixation and permeabilization reagents. Here, we outline protocols for the staining of lipids in different cellular compartments. Please read the entire protocol before beginning.**

### **Procedure, general lipid staining:**

1. Fix cells by adding 4% paraformaldehyde to cells in media for 20 minutes at room temperature.
2. Wash three times with TBS.
3. Permeabilize the cells with 0.5% Saponin or 0.01% Digitonin at RT for 15 minutes.
4. Wash three times with TBS.
5. Block with 10% Goat Serum in TBS either overnight at 4 °C or 30 minutes at 37 °C.
6. Add the anti-lipid or anti-PIP antibody diluted in 10% Goat Serum in TBS to the concentration suggested on the technical data sheet. Incubate for 60 minutes at 37 °C.
7. Remove the primary antibody solution and wash for 5-10 minutes in TBS-Goat Serum 1%. Repeat twice.
8. Dilute the secondary antibody in 10% Goat Serum in TBS to the concentration suggested on the technical data sheet. Add the secondary antibody solution and incubate for 45 minutes.
  - For fluorescently conjugated secondary antibodies, this step and all remaining steps until mounting should be performed covered or in the dark.
9. Remove the secondary antibody solution and wash for 5-10 minutes in TBS. Repeat twice.
10. Remove excess wash buffer by blotting the edge of the coverslip.
  - DAPI or alternative nucleic acid stains may also be applied at this time if they are not present in the mounting media. Excess nucleic acid stain should also be removed prior to mounting.
11. Add the mounting media to a microscope slide and mount and seal the coverslip. Repeat this for all stained samples and store at 4 °C in the dark until imaging.

### **Procedure, plasma membrane staining:**

1. Remove culture media from cells and fix with 4% formaldehyde (FA) and 0.2% glutaraldehyde (GA) in PBS for 15 minutes at room temperature (RT).
  - Fixation may also be performed at 4 °C for 3 hours.**Steps 2-8 should be performed on ice or at 4 °C using pre-chilled buffers and solutions.**
2. Remove the fixation solution and rinse three times with 50 mM NH<sub>4</sub>Cl in PBS.
3. Block and permeabilize the cells with Buffer 1 (see recipe below) supplemented with 5% normal goat serum, 0.5% saponin, and 50 mM NH<sub>4</sub>Cl for 45 minutes.
4. Dilute the primary antibody in Buffer 2 (see recipe below) to the concentration suggested on the technical data sheet. Remove the block and permeabilization solution and add the primary antibody solution. Incubate for 60 minutes.
5. Remove the primary antibody solution and wash for 5-10 minutes in Buffer 1. Repeat once.
6. Dilute the secondary antibody in Buffer 2 to the concentration suggested on the technical data sheet. Add the secondary antibody solution and incubate for 45 minutes.
  - For fluorescently conjugated secondary antibodies, this step and all remaining steps until mounting should be performed covered or in the dark.
7. Rinse the slides or coverslips with Buffer 1 four times.
8. Post fix cells in 2% FA in PBS for 10 minutes.
9. Following post-fixation, bring the cells to RT. All remaining steps may be performed at RT.
10. Rinse cells three times with 50 mM NH<sub>4</sub>Cl in PBS, then once with MQH<sub>2</sub>O or dH<sub>2</sub>O.
11. Remove excess wash buffer by blotting the edge of the coverslip.
  - DAPI or alternative nucleic acid stains may also be applied at this time if they are not present in the mounting media. Excess nucleic acid stain should also be removed prior to mounting.

# Echelon Biosciences Inc.

---

12. Add the mounting media to a microscope slide and mount and seal the coverslip. Repeat this for all stained samples and store at 4 °C in the dark until imaging.

Buffer 1: (20 mM PIPES, pH 6.8, 137 mM NaCl, 2.7 mM KCl)

Buffer 2: (20 mM PIPES, pH 6.8, 137 mM NaCl, 2.7 mM KCl) + 5% normal goat serum, 0.1% saponin

## **Procedure. Golgi apparatus staining:**

**All steps should be performed at room temperature.**

1. Remove culture media from cells and fix with 2% FA in PBS for 15 minutes.
2. Remove the fixation solution and rinse three times with 50 mM NH<sub>4</sub>Cl in PBS.
3. Permeabilize the cells with Buffer 1 supplemented with 20 µM digitonin for 5 minutes.
4. Rinse three times with Buffer 1 and block the cells in Buffer 1 supplemented with 5% normal goat serum and 50 mM NH<sub>4</sub>Cl for 45 minutes.
5. Dilute the primary antibody in Buffer 2 to the concentration suggested on the technical data sheet. Remove the block and permeabilization solution and add the primary antibody solution. Incubate for 60 minutes.
6. Remove the primary antibody solution and wash for 5-10 minutes twice in Buffer 1.
7. Dilute the secondary antibody in Buffer 2 to the concentration suggested on the technical data sheet. Add the secondary antibody solution and incubate for 45 minutes.
  - For fluorescently conjugated secondary antibodies, this step and all remaining steps until mounting should be performed covered or in the dark.
8. Rinse the slides or coverslips with Buffer 1 four times.
9. Post fix cells in 2% FA in PBS for 10 minutes.
10. Rinse cells three times with 50 mM NH<sub>4</sub>Cl in PBS, then once with MQH<sub>2</sub>O or dH<sub>2</sub>O.
11. Remove excess wash buffer and dry cells.
  - DAPI or alternative nucleic acid stains may also be applied at this time if they are not present in the mounting media. Excess nucleic acid stain should also be removed prior to mounting.
12. Add the mounting media to a microscope slide and mount and seal the coverslip. Repeat this for all stained samples and store at 4 °C in the dark until imaging.

Buffer 1: (20 mM PIPES, pH 6.8, 137 mM NaCl, 2.7 mM KCl)

Buffer 2: Buffer 1 + 5% normal goat serum.

## **Procedure. nuclear membrane staining:**

**All steps should be performed at room temperature.**

For staining of the nuclear membrane, please refer to and follow the protocol described for 'Golgi apparatus staining' and substitute 0.2% Triton X-100 for digitonin.

## **Additional Notes**

1. These protocols are recommended as starting points and are based on published data for lipid and phosphatidylinositol phosphate immunostaining. Users are encouraged to adjust the protocols as needed based on the provided references.
2. Variability in staining patterns may be observed across different cell types owing to distribution of specific lipids and cellular lipid composition.
3. The stringency of the fixation conditions can be adjusted by increasing the crosslinker concentration, i.e. 2% FA --> 4% FA --> 4% FA + 0.2% GA, as well as the temperature.
4. Most common crosslinking reagents are unlikely to react with lipids directly and will instead react mainly with proteins. As such, performing staining at 4 °C vs room temperature may affect residual lipid mobility after an initial fixation in some membranes.
5. Post-fixation of cells after antibody staining may also reduce residual lipid mobility.
6. In general, we do not recommend the use of acetone or methanol as cellular fixatives as they are organic solvents and can solubilize and extract lipids.
7. If the use of acetone or methanol is required to visualize a protein of interest during co-immunostaining for a specific lipid, then we encourage the user to compare the staining pattern for a given lipid with an alternative fixation procedure.
8. If using the 'general lipid staining' protocol, the choice of permeabilization reagent should be made with the target



# Echelon Biosciences Inc.

cellular compartment in mind: digitonin and saponin are selective for cholesterol and therefore require it for effective permeabilization. Triton X-100 and Tween20 will permeabilize all membranes but may also completely solubilize some membranes such as the Golgi and ER.

9. At sufficiently high concentrations, Saponin, Triton X-100, and Tween20 may lyse cells or extract lipids. We do not recommend exceeding 0.5% Saponin, 0.2% Triton X-100, or 0.1% Tween20.

## Protocol References

1. Hammond, G.R., et al., PI4P and PI(4,5)P2 are essential but independent lipid determinants of membrane identity. *Science*, 2012. 337(6095): p. 727-30.
2. Hammond, G.R., G. Schiavo, and R.F. Irvine, Immunocytochemical techniques reveal multiple, distinct cellular pools of PtdIns4P and PtdIns(4,5)P(2). *Biochem J*, 2009. 422(1): p. 23-35.
3. Hammond, G.R., et al., Elimination of plasma membrane phosphatidylinositol (4,5)-bisphosphate is required for exocytosis from mast cells. *J Cell Sci*, 2006. 119(Pt 10): p. 2084-94.

## Antibody References

1. Qin L, Zhou Z, Li Q, Zhai C, Liu L, Quilichini TD, et al. Specific Recruitment of Phosphoinositide Species to the Plant-Pathogen Interfacial Membrane Underlies Arabidopsis Susceptibility to Fungal Infection. *The Plant Cell*. 2020
2. Pan P-Y, Sheehan P, Wang Q, Zhu X, Zhang Y, Choi I, et al. Synj1 Haploinsufficiency Causes Dopamine Neuron Vulnerability and alpha-synuclein Accumulation in Mice. *Human Molecular Genetics*. 2020.
3. Palušová V, Renzová T, Verlande A, Vaclová T, Medková M, Cetlová L, et al. Dual Targeting of BRAF and mTOR Signaling in Melanoma Cells with Pyridinyl Imidazole Compounds. *Cancers*. 2020;12(6):1516.
4. Fujii Y, Murata-Kamiya N, Hatakeyama M. Helicobacter pylori CagA oncoprotein interacts with SHIP2 to increase its delivery into gastric epithelial cells. *Cancer Science*. 2020;111(5):1596-606.
5. Fáberová V, Kalasová I, Krausová A, Hozák P. Super-Resolution Localisation of Nuclear PI(4)P and Identification of Its Interacting Proteome. *Cells*. 2020;9(5):1191.
6. Amiar S, Katris NJ, Berry L, Dass S, Duley S, Arnold C-S, et al. Division and Adaptation to Host Environment of Apicomplexan Parasites Depend on Apicoplast Lipid Metabolic Plasticity and Host Organelle Remodeling. *Cell Reports*. 2020;30(11):3778-92.e9.
7. Wang H, Tai AW. Nir2 Is an Effector of VAPs Necessary for Efficient Hepatitis C Virus Replication and Phosphatidylinositol 4-Phosphate Enrichment at the Viral Replication Organelle. *Journal of virology*. 2019;93(22):e00742-19.
8. Wang H, Loerke D, Bruns C, Müller R, Koch PA, Puchkov D, et al. Phosphatidylinositol 3,4-bisphosphate synthesis and turnover are spatially segregated in the endocytic pathway. *Journal of Biological Chemistry*. 2019.
9. Ramos AR, Ghosh S, Suhel T, Chevalier C, Obeng EO, Fafilek B, et al. Phosphoinositide 5-phosphatases SKIP and SHIP2 in ruffles, the endoplasmic reticulum and the nucleus: an update. *Advances in Biological Regulation*. 2019:100660.
10. Park MK, Yao Y, Xia W, Setijono SR, Kim JH, Vila IK, et al. PTEN self-regulates through USP11 via the PI3K-FOXO pathway to stabilize tumor suppression. *Nature Communications*. 2019;10(1):636.
11. Miao G, Zhang Y, Chen D, Zhang H. The ER-Localized Transmembrane Protein TMEM39A/SUSR2 Regulates Autophagy by Controlling the Trafficking of the PtdIns(4)P Phosphatase SAC1. *Molecular Cell*. 2019.
12. Mahajan D, Tie HC, Chen B, Lu L. Dopey1-Mon2 complex binds to dual-lipids and recruits kinesin-1 for membrane trafficking. *Nature Communications*. 2019;10(1):3218.
13. Li Z, Guo D, Qin Y, Chen M. PI4KB on Inclusion Bodies Formed by ER Membrane Remodeling Facilitates Replication of Human Parainfluenza Virus Type 3. *Cell Reports*. 2019;29(8):2229-42.e4. doi: <https://doi.org/10.1016/j.celrep.2019.10.052>.
14. Koike S, Jahn R. SNAREs define targeting specificity of trafficking vesicles by combinatorial interaction with tethering factors. *Nature Communications*. 2019;10(1):1608.
15. Judith D, Jefferies HBJ, Boeing S, Frith D, Snijders AP, Tooze SA. ATG9A shapes the forming autophagosome through Arfaptin 2 and phosphatidylinositol 4-kinase III $\beta$ . *The Journal of cell biology*. 2019;218(5):1634-52. Epub 03/27.
16. Horova V, Lyoo H, Rózycki B, Chalupska D, Smola M, Humpolickova J, et al. Convergent evolution in the mechanisms of ACBD3 recruitment to picornavirus replication sites. *PLOS Pathogens*. 2019;15(8):e1007962.
17. Feng Z, Xu K, Kovalev N, Nagy PD. Recruitment of Vps34 PI3K and enrichment of PI3P phosphoinositide in the viral replication compartment is crucial for replication of a positive-strand RNA virus. *PLOS Pathogens*. 2019;15(1):e1007530.
18. 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- $\kappa$ B via phosphoinositide remodeling of endosomes. *Nature Communications*. 2019;10(1):2247.
19. Du X, Zhou L, Aw YC, Mak HY, Xu Y, Rae J, et al. ORP5 localizes to ER-lipid droplet contacts and regulates the level of PI(4)P on lipid droplets. *The Journal of Cell Biology*. 2019;219(1).
20. Dong J, Du X, Wang H, Wang J, Lu C, Chen X, et al. Allosteric enhancement of ORP1-mediated cholesterol transport by PI(4,5)P2/PI(3,4)P2. *Nature Communications*. 2019;10(1):829.



# Echelon Biosciences Inc.

---

21. Di Pardo A, Pepe G, Castaldo S, Marracino F, Capocci L, Amico E, et al. Stimulation of Sphingosine Kinase 1 (SPHK1) Is Beneficial in a Huntington's Disease Pre-clinical Model. *Frontiers in Molecular Neuroscience*. 2019;12(100).
22. Choi S, Chen M, Cryns VL, Anderson RA. A nuclear phosphoinositide kinase complex regulates p53. *Nature Cell Biology*. 2019;21:462-75.
23. Wang H, Ma Q, Qi Y, Dong J, Du X, Rae J, et al. ORP2 Delivers Cholesterol to the Plasma Membrane in Exchange for Phosphatidylinositol 4, 5-Bisphosphate (PI(4,5)P2). *Molecular Cell*. 2018.
24. 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.
25. Sbrissa D, Semaan L, Govindarajan B, Li Y, Caruthers NJ, Stemmer PM, et al. A novel cross-talk between CXCR4 and PI4KIII $\alpha$  in prostate cancer cells. *Oncogene*. 2018.

