

Cell lines & organisms analyzed with Propargyl-choline (PCho) and 1-Azidoethyl-choline (AECho)

Cell line / Organism	Final PCho concentration	Final AECho concentration
NIH3T3 cells	10 μ M – 5 mM ^[1] , 100 μ M ^[2]	250 – 500 μ M ^[2]
HEK293T cells	100 μ M ^[1]	
Jurkat cells	300 – 500 μ M ^[3] , 400 μ M ^[7]	300 – 500 μ M ^[3]
HeLa cells	200 μ M ^[6] , 2 mM ^[8]	75 μ M ^[3]
Vero cells	12.5 μ M – 1.6 mM ^[4]	12.5 μ M – 1.6 mM ^[4] 400 μ M ^[5]
PC12 cells, clone 251	400 μ M ^[7]	
HepG2 cells	400 μ M ^[7]	
BT-474 cells	2 mM ^[8]	
MDA-MB-231 cells	2 mM ^[8]	
5637-sensitive cells	2 mM ^[8]	
5637-resistant cells	2 mM ^[8]	
Mouse	1 M (50 μ mol) ^[1]	

Selected References:

- [1] Jao *et al.* (2009) Metabolic labeling and direct imaging of choline phospholipids in vivo. *Proc. Natl. Acad. Sci. USA* **106(36)**: 15332.
- [2] Jao *et al.* (2015) Biosynthetic Labeling and Two-Color Imaging of Phospholipids in Cells. *Chem. Bio. Chem.* **16**:472.
- [3] Caishun *et al.* (2014) Practical Labeling Methodology for Choline-Derived Lipids and Applications in Live Cell Fluorescence Imaging. *Photochemistry and Photobiology* **90**: 686.
- [4] Huang *et al.* (2013) Enveloped Virus Labeling via Both Intrinsic Biosynthesis and Metabolic Incorporation of Phospholipids in Host Cells. *Anal. Chem.* **85**: 5236.
- [5] Zhao *et al.* (2016) Surface labeling of enveloped virus with polymeric imidazole ligand-capped quantum dots via the metabolic incorporation of phospholipids into host cells. *J. Mater. Chem. B* **4**: 2421.
- [6] Zhang *et al.* (2016) Positive-strand RNA viruses stimulate host phosphatidylcholine synthesis at viral replication sites. *Proc. Natl. Acad. Sci. USA* **113(8)**: E1064.
- [7] Merklinger *et al.* (2016) No Evidence for Spontaneous Lipid Transfer at ER-PM Membrane Contact Sites. *J. Membrane. Biol.* **249**: 41.
- [8] Luo *et al.* (2016) Rapid assessment of drug resistance of cancer cells to gefitinib and carboplatin using optical imaging. *Anal. Biochem.* **504**: 50.
- [9] Luo *et al.* (2013) High-Resolution Optical Molecular Imaging of Changes in Choline Metabolism in Oral Neoplasia. *Translational Oncology* **6**: 33.