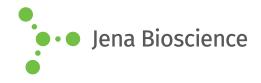
DATA SHEET

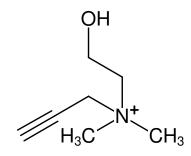




Propargyl-choline

Bromo salt

Cat. No.	Amount
CLK-066	1 mg



Structural formula of Propargyl-choline

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient

 $temperature\ possible.$

Shelf Life: 24 months after date of delivery

Molecular Formula: C₇H₁₄NO (cation)
Molecular Weight: 128.19 g/mol (cation)

Purity: H NMR conforms to structure

Form: solid or clear oil **Color:** white to off-white

Solubility: water, 1x PBS buffer

Applications:

Choline-containing phospholipid synthesis monitoring in cell culture and whole organisms^[1-4]

Related Products:

Copper (II)-Sulphate (CuSO₄), #CLK-MI004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-MI005 1-Azidoethyl-choline, #CLK-065

Selected References:

[1] Jao *et al.* (2009) Metabolic Labeling and direct imaging of choline phospholipids in vivo. *PNAS* **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] Zhao *et al.* (2016) Surface labeling of enveloped virus with polymeric imidazole-capped quantum dots via the metabolic incorporation of phospholipids into host cells. *J. Mater. Chem. B* **4**:2421.

Di Guilmi *et al.* (2017) Specific and spatial labeling of choline-containing teichoic acids in *Streptococcus pneumoniae* by click chemistry *Chem. Commun.* 10.1039/c7cc05646j.