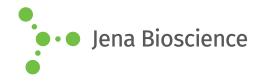
# **DATA SHEET**





## 1-Azidoethyl-choline

Iodo salt

Cat. No.	Amount
CLK-065	1 mg

Structural formula of 1-Azidoethyl-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:** C<sub>6</sub>H<sub>15</sub>N<sub>4</sub>O (cation) **Molecular Weight:** 159.21 g/mol (cation)

Exact Mass: 159.12 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<sup>[1-4]</sup>

## **Related Products:**

Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-MI004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-MI005 Propargyl-choline, #CLK-066

#### Selected References:

[1] Jao et al. (2015) Biosynthetic Labeling and Two-Color Imaging of Phospholipids in Cells. Chem. Bio. Chem. **16**:472.

[2] Caishun et al. (2014) Practical Labeling Methodology for Choline-derived Lipids and Applications in Live Cell Fluorescence Imaging. *Photochemistry and Photobiology* **90**:686.

[3] Huang *et al.* (2013) Enveloped Virus Labeling via Both Intrinsic Biosynthesis and Metabolic Incorporation of Phospholipids in Host Cells. *Anal. Chem.* **85**:5236.

[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.