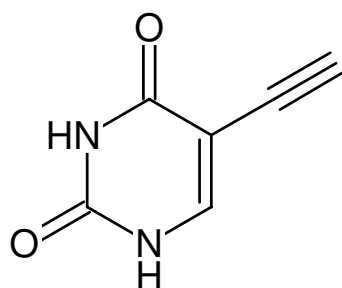


**5-Ethynyluracil**

Cat. No.	Amount
CLK-089-10	10 mg



Structural formula of 5-Ethynyluracil

Selected References:

Zajaczkowski *et al.* (2018) Bioorthogonal Metabolic Labeling of Nascent RNA in Neurons Improves the Sensitivity of Transcriptome-Wide Profiling. *ACS Chem. Neurosci.* **9** (7):1858.

For research use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₆H₄N₂O₂

Molecular Weight: 136.11 g/mol

Exact Mass: 136.03 g/mol

CAS#: 59989-18-3

Purity: ≥ 97 % (HPLC)

Form: solid

Color: white to off-white

Solubility: DMSO, MeOH