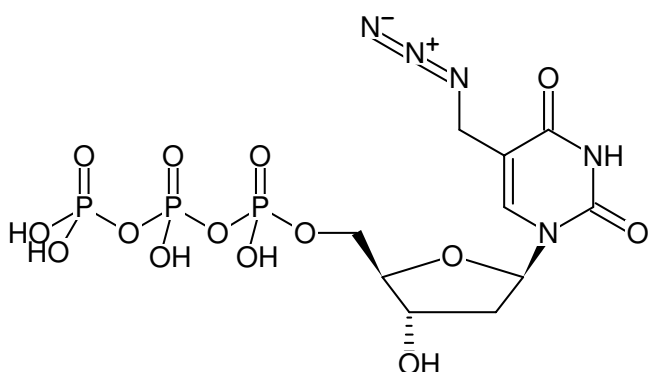


**5-Azidomethyl-dUTP**

5-AmdUTP

5-Azidomethyl-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-084S	20 µl (10 mM)
CLK-084L	5 x 20 µl (10 mM)



Structural formula of 5-Azidomethyl-dUTP

For research use only!**Shipping:** shipped on blue ice**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:** C₁₀H₁₆N₅O₁₄P₃ (free acid)**Molecular Weight:** 523.18 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 267 nm; ε 9.65 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)