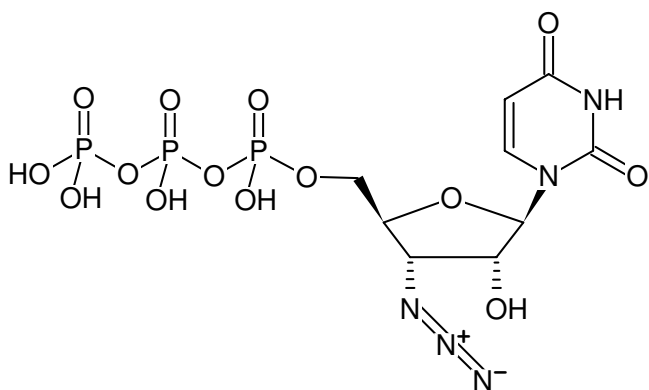


**3'-Azido-3'-dUTP**

3'-Azido-3'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-992	100 µl (10 mM)



Structural formula of 3'-Azido-3'-dUTP

For research use only!**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: 12 months after date of delivery**Molecular Formula:** C₉H₁₄N₅O₁₄P₃ (free acid)**Molecular Weight:** 509.15 g/mol (free acid)**Exact Mass:** 508.98 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 262 nm, ε 9.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)