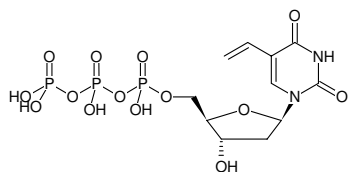


**5-Vinyl-dUTP**

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

Cat. No.	Amount
CLK-068	50 µl (10 mM)



Structural formula of 5-Vinyl-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:** C₁₁H₁₇N₂O₁₄P₃ (free acid)**Molecular Weight:** 494.18 g/mol (free acid)**Exact Mass:** 493.99 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 292 nm, ε 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)