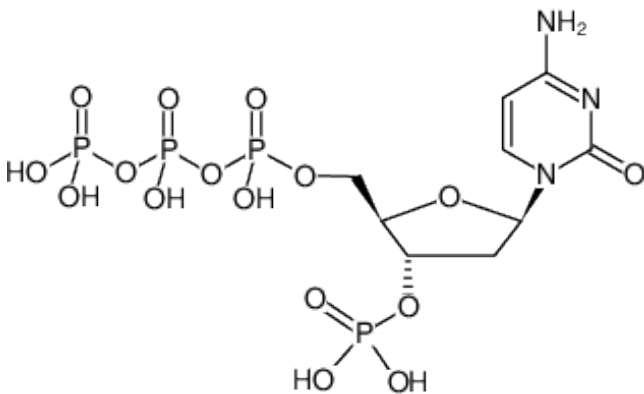


**3'-Phosphate-dCTP**

3'-Phosphate-2'-deoxycytidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1222S	200 µl (10 mM)
NU-1222L	5 x 200 µl (10 mM)



Structural formula of 3'-Phosphate-dCTP

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: 12 months after date of delivery**Molecular Formula:** C₉H₁₇N₃O₁₆P₄ (free acid)**Molecular Weight:** 547.14 g/mol (free acid)**Exact Mass:** 546.96 g/mol (free acid)**Purity:** ≥ 98 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)