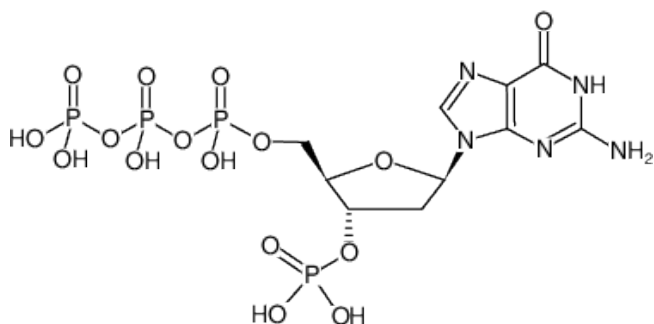




3'-Phosphate-dGTP

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

Cat. No.	Amount
NU-1223S	200 µl (10 mM)
NU-1223L	5 x 200 µl (10 mM)



Structural formula of 3'-Phosphate-dGTP

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: C₁₀H₁₇N₅O₁₆P₄ (free acid)

Molecular Weight: 587.16 g/mol (free acid)

Exact Mass: 586.96 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} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)