

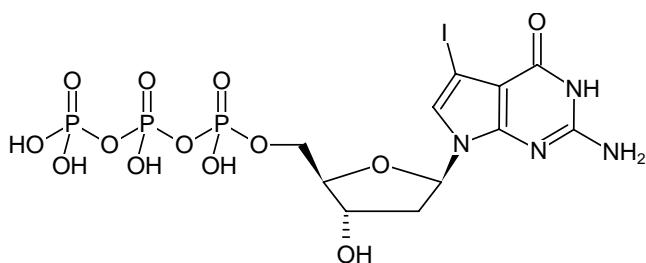


7-Deaza-7-iodo-dGTP

(7-Deaza-7-I-dGTP)

7-Deaza-7-iodo-2'-deoxyguanosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1182S	50 µl (10 mM)
NU-1182L	5 x 50 µl (10 mM)



Structural formula of 7-Deaza-7-iodo-dGTP

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: C₁₁H₁₆N₄O₁₃P₃I (free acid)

Molecular Weight: 632.09 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 11.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)