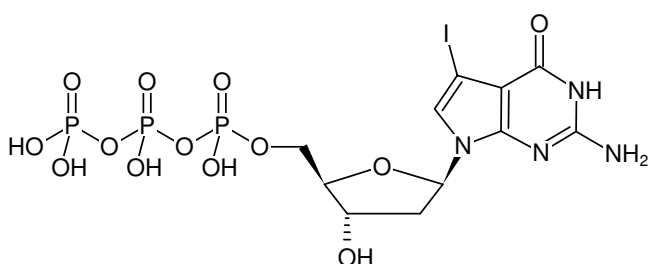


**7-Deaza-7-iodo-dGTP**

(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 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:** C₁₁H₁₆N₄O₁₃P₃I (free acid)**Molecular Weight:** 632.09 g/mol (free acid)**Exact Mass:** 631.90 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} 267 nm, ε 11.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)