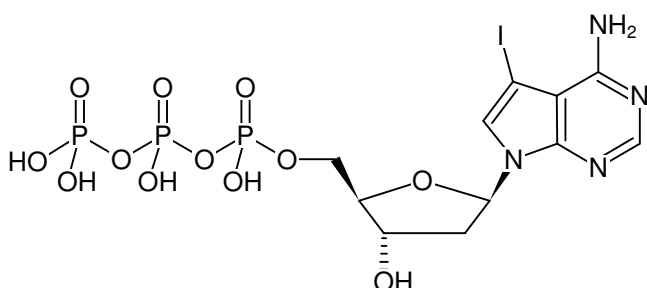


**7-Deaza-7-iodo-dATP**

(7-Deaza-7-I-dATP)

7-Deaza-7-iodo-2'-deoxyadenosine-5'-triphosphate, Sodium salt

| Cat. No. | Amount |
|----------|-------------------|
| NU-1181S | 50 µl (10 mM) |
| NU-1181L | 5 x 50 µl (10 mM) |



Structural formula of 7-Deaza-7-iodo-dATP

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:** C₁₁H₁₆N₄O₁₂P₃I (free acid)**Molecular Weight:** 616.09 g/mol (free acid)**Exact Mass:** 615.90 g/mol (free acid)**CAS#:** 214833-26-8**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 283 nm, ε 8.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)