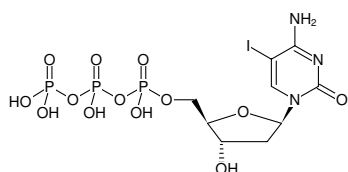


**5-Iodo-dCTP**

(5I-dCTP)

5-Iodo-2'-deoxycytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-128S	50 µl (10 mM)
NU-128L	5 x 50 µl (10 mM)



Structural formula of 5-Iodo-dCTP

For research use only!**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:** C₉H₁₅N₃O₁₃P₃I (free acid)**Molecular Weight:** 593.05 g/mol (free acid)**Exact Mass:** 592.89 g/mol (free acid)**CAS#:** 31747-59-8**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} 293 nm, ε 5.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)