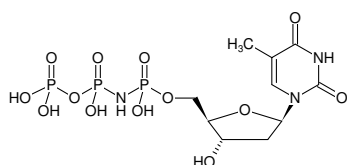


**dTpNHpp**

(dTMPNPP)

2'-Deoxythymidine-5'-[(α,β)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-907-1	1 mg



Structural formula of dTpNHpp

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:** C₁₀H₁₈N₃O₁₃P₃ (free acid)**Molecular Weight:** 481.18 g/mol (free acid)**Exact Mass:** 481.01 g/mol (free acid)**CAS#:** 141171-14-4 (free acid), 183584-80-7 (Tetrasodium salt)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 267 nm, ϵ 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Koag *et al.* (2019) Mutagenic Replication of the Major Oxidative Adenine Lesion 7,8-Dihydro-8-oxoadenine by Human DNA Polymerases. *J. Am. Chem. Soc.* **141** (11):4584.

Ma *et al.* (2008) New thymidine triphosphate analogue inhibitors of human immunodeficiency virus-1 reverse transcriptase. *J. Med. Chem.* **35** (11):1938.