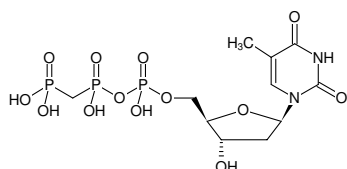


**dTppCp**

(dTMPPCP)

2'-Deoxythymidine-5'-[β,γ]-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-262	5 mg



Structural formula of dTppCp

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:** C₁₁H₁₉N₂O₁₃P₃ (free acid)**Molecular Weight:** 480.20 g/mol (free acid)**Exact Mass:** 480.01 g/mol (free acid)**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)