

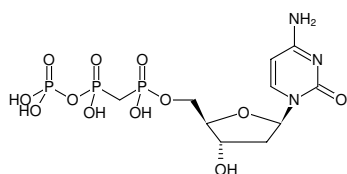
**dCpCpp**

(dCMP CPP)

dCpCpp (dCMP CPP)

2'-Deoxycytidine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-437S	50 $\mu$ l (10 mM)
NU-437L	5 x 50 $\mu$ l (10 mM)



Structural formula of dCpCpp

**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:** C<sub>10</sub>H<sub>18</sub>N<sub>3</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 465.18 g/mol (free acid)**Exact Mass:** 465.01 g/mol (free acid)**CAS#:** 941569-33-1**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  271 nm,  $\epsilon$  9.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**

Vaidyanathan *et al.* (2013) Insights into the Conformation of Aminofluorene-Deoxyguanine Adduct in a DNA Polymerase Active Site *J. Biol. Chem.* **288** (32):23573–23585.

Liang *et al.* (2008) Alpha,beta-methylene-2'-deoxynucleoside 5'-triphosphates as noncleavable substrates for DNA polymerases: isolation, characterization, and stability studies of novel 2'-deoxycyclonucleosides, 3,5'-cyclo-dG, and 2,5'-cyclo-dT. *J. Med. Chem.* **51** (20):6460-70.