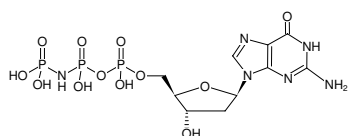


**dGppNHp**

(dGMPPNP)

2'-Deoxyguanosine-5'-[( $\beta,\gamma$ )-imido]triphosphate, Lithium salt

Cat. No.	Amount
NU-258	5 mg



Structural formula of dGppNHp

**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:** 6 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>17</sub>N<sub>6</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 506.20 g/mol (free acid)**Exact Mass:** 506.01 g/mol (free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\max}$  252 nm,  $\epsilon$  13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)