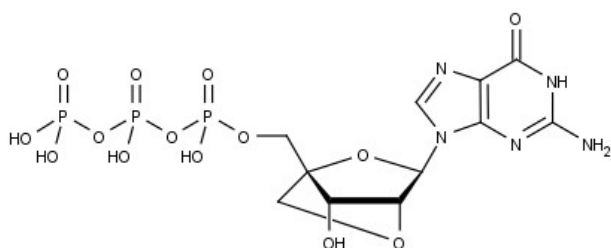


**LNA-GTP**

LNA-guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-983	10 µl (100 mM)



Structural formula of LNA-GTP

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:** C₁₁H₁₆N₅O₁₄P₃ (free acid)**Molecular Weight:** 535.19 g/mol (free acid)**Exact Mass:** 534.99 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Related Products:**

HighFidelity LNA PCR Labeling Kit, #APP-101-LNA

Selected References:Veedu *et al.* (2010) Polymerase directed incorporation studies of LNA-G nucleoside 5'-triphosphate and primer extension involving all four LNA nucleotides. *New J. Chem.* **34**:877.