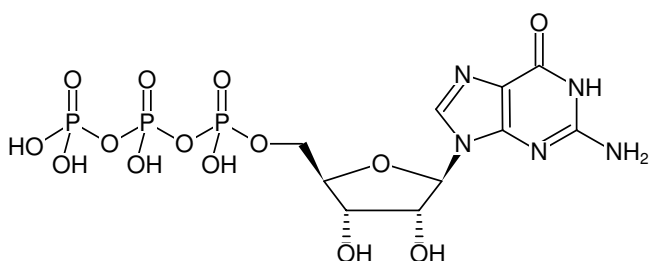


**GTP - Solid**

Guanosine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1047-200	200 mg
NU-1047-1G	1 g
NU-1047-10G	10 g
NU-1047-100G	100 g

Quality Control Specifications: in vitro transcription (T7 RNA polymerase): visible RNA fragments after 5 min incubation, Proteases: not detectable



Structural formula of GTP - Solid

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

Molecular Formula: C₁₀H₁₆N₅O₁₄P₃ (free acid)

Molecular Weight: 523.18 g/mol (free acid)

Exact Mass: 522.99 g/mol (free acid)

CAS#: 36051-31-7

Purity: ≥ 90 % (HPLC)

Form: Powder

Color: White to slightly yellow

Solubility: Soluble in water. Clear solution.

Spectroscopic Properties: λ_{max} 252 nm, ε 14.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)