



GP₃G (Unmethylated Cap Analog) - Solid

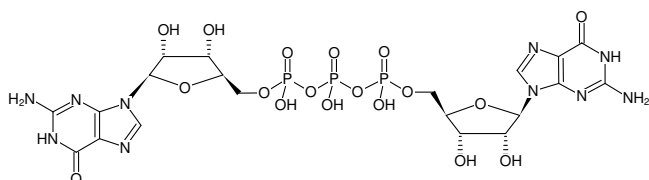
G(5')ppp(5')G

P¹-(Guanosyl) P³-(5'-(guanosyl))triphosphate, Sodium salt

Cat. No.	Amount
NU-854-1	1 mg
NU-854-5	5 mg

Selected References:

Liu *et al.* (2002) The scavenger mRNA decapping enzyme DcpS is a member of the HIT family of pyrophosphatases. *EMBO J.* **21** (17):397.



Structural formula of GP₃G (Unmethylated Cap Analog) - 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 after date of delivery

Molecular Formula: C₂₀H₂₇N₁₀O₁₈P₃ (free acid)

Molecular Weight: 788.41 g/mol (free acid)

Exact Mass: 788.07 g/mol (free acid)

CAS#: 6674-45-9 (acid), 1314754-91-0 (Trisodium salt)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 251 nm, ε 23.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)