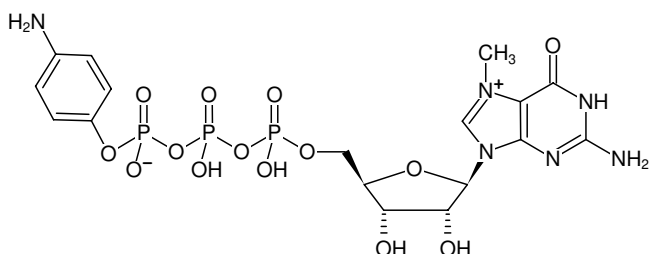




γ -Aminophenyl-m⁷GTP

7-Methylguanosine-5'-[γ -(4-aminophenyl)]triphosphate, Sodium salt

| Cat. No. | Amount |
|----------|--------|
| NU-870-1 | 1 mg |



Structural formula of γ -Aminophenyl-m⁷GTP

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: C₁₇H₂₃N₆O₁₄P₃ (free acid)

Molecular Weight: 628.32 g/mol (free acid)

Exact Mass: 628.05 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{\max} 260 nm, ϵ 10.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)