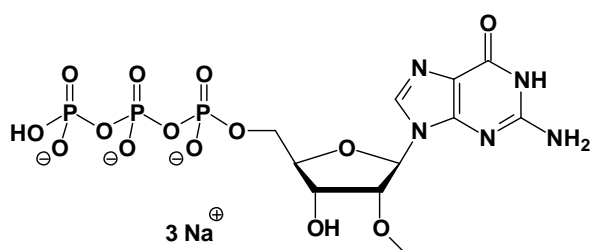


2'-O-Methyl-GTP

2'-O-Methylguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1127S	150 Units
NU-1127L	750 Units



Cat. No.: NU-1127

Molecular Formula: C₁₁H₁₅N₅O₁₄P₃ (Anion)

Molecular Weight: 534.18 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{\max} 252 nm;
 ϵ 13700

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Fresno *et al.* (1980) Inhibitory effects of 'cap' analogues on globin mRNA and encephalomyocarditis RNA translation in a reticulocyte cell-free system. *Eur J Biochem.* **103(1)**:125.