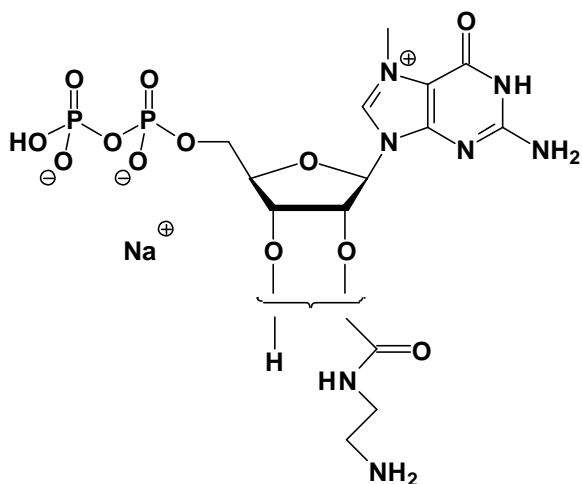


EDA-m⁷-GDP

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate,
Sodium salt

Cat. No.	Amount
NU-827S	30 Units
NU-827L	150 Units



Cat. No.: NU-827

Molecular Formula: C₁₄H₂₂N₇O₁₂P₂ (Anion)

Molecular Weight: 542.33 (Anion)

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

Spectroscopic properties: λ_{max} 258/280 nm;
ε 9800/8000

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 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution