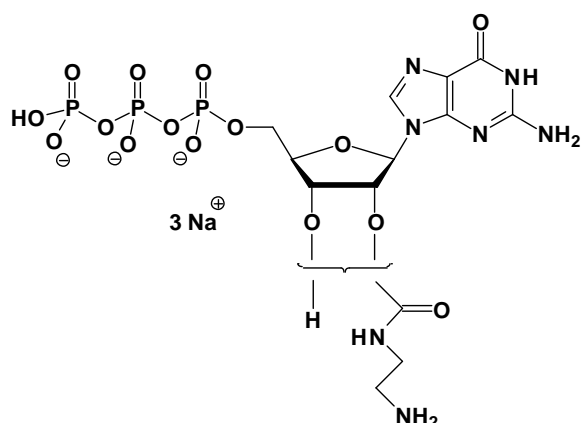


EDA-GTP

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-820S	30 Units
NU-820L	150 Units



Cat. No.: NU-820

Molecular Formula: C₁₃H₁₉N₇O₁₅P₃ (Anion)

Molecular Weight: 606.25 (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 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution