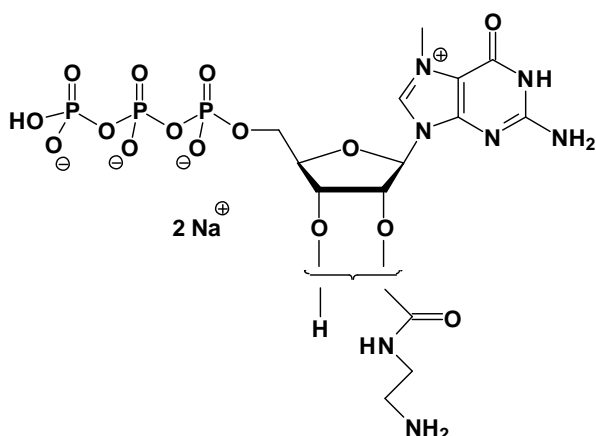


## EDA-m<sup>7</sup>GTP

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

Cat. No.	Amount
NU-824S	30 Units
NU-824L	150 Units



**Cat. No.:** NU-824

**Molecular Formula:** C<sub>14</sub>H<sub>22</sub>N<sub>7</sub>O<sub>15</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 621.31 (Anion)

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

**Spectroscopic properties:** λ<sub>max</sub> 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