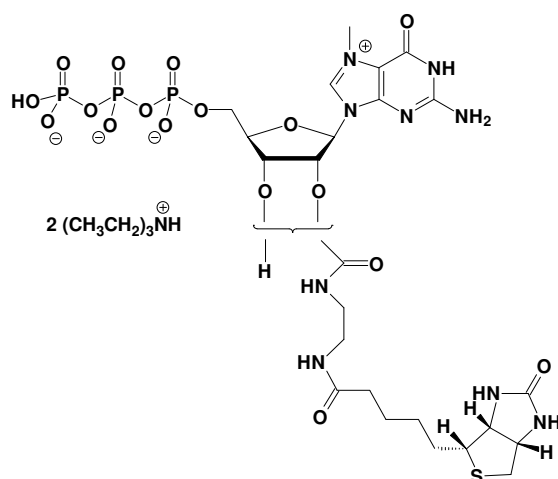


EDA-m⁷GTP - Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-N7-methyl-guanosine-5'-triphosphate - Biotin,
Triethylammonium salt

Cat. No.	Amount
NU-824-BIO	50 μ l / 5 mM



Cat. No.: NU-824-BIO

Molecular Formula: C₂₄H₃₆N₉O₁₇P₃S (Anion)

Molecular Weight: 847.58 (Anion)

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

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!