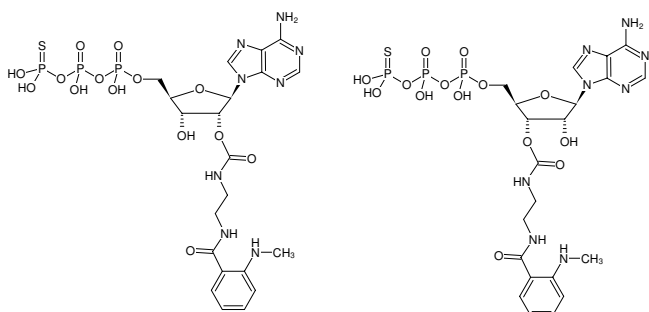


**EDA-ATPyS-MANT**

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-1609-MNT	50 µl (1 mM)



Structural formula of EDA-ATPyS-MANT

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₂₁H₂₉N₈O₁₄P₃S (free acid)

Molecular Weight: 742.48 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 335 nm, λ_{em} 440 nm, ε 2.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)