



EDA-ATPyS-ATTO-612Q

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with ATTO 612Q, Triethylammonium salt

Cat. No.	Amount
NU-1609-612Q	50 µl (1 mM)

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: C₁₃H₂₂N₇O₁₃P₃S - ATTO 612Q (free acid)

Molecular Weight: 1281.33 g/mol (free acid)

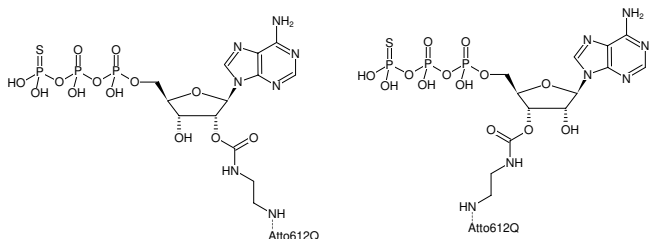
Purity: ≥ 90 % (HPLC)

Form: solution in water

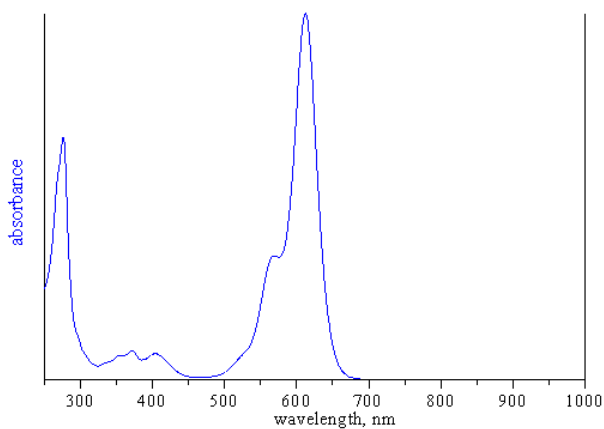
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 615 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ATPyS-ATTO-612Q



absorption spectrum of ATTO 612Q