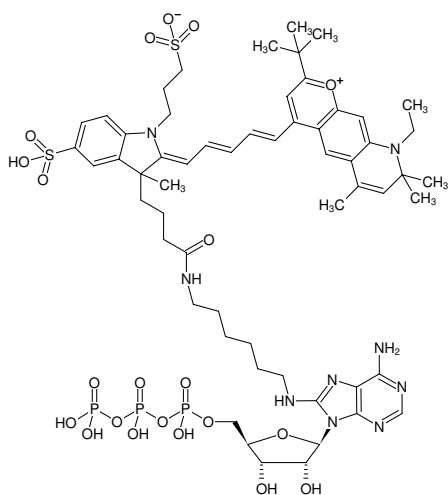




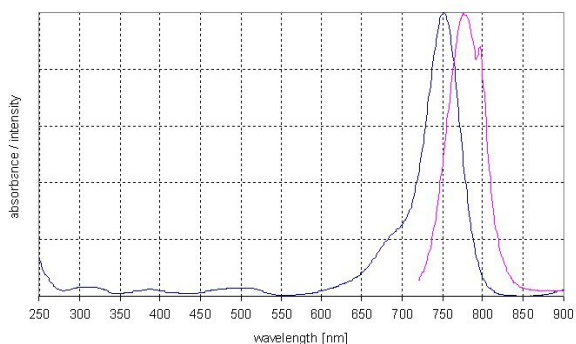
8-(6-Aminoethyl)-amino-ATP-DY-751

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, labeled with DY 751, Triethylammonium salt

Cat. No.	Amount
NU-807-751	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-ATP-DY-751



excitation and emission spectrum of DY 751

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₈H₈₀N₉O₂₁P₃S₂ (free acid)

Molecular Weight: 1395.41 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 751 nm, λ_{em} 779 nm, ε 270.0 L mol⁻¹ cm⁻¹ (Ethanol)