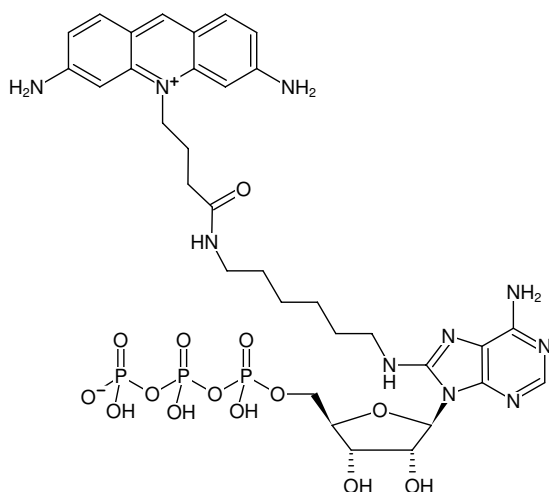




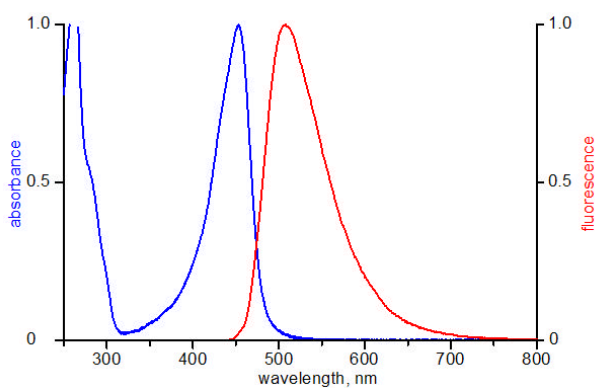
8-(6-Aminoethyl)-amino-ATP-ATTO-465

8-(6-Aminoethyl)-amino-adenosine-5'-triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-807-465	120 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-ATP-ATTO-465



excitation and emission spectrum of ATTO 465

For general laboratory use.

Shipping: shipped on gel packs

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: C₃₃H₄₅N₁₀O₁₄P₃ (free acid)

Molecular Weight: 898.69 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 453 nm, λ_{em} 506 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)