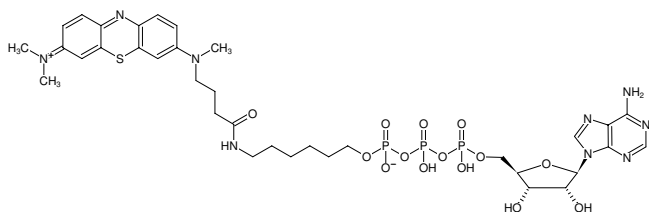




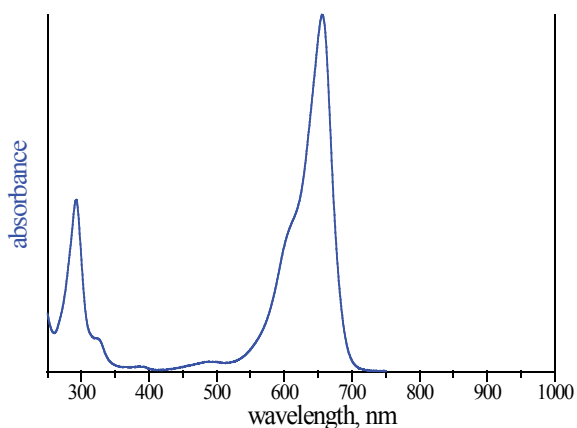
γ -(6-Aminoethyl)-ATP-ATTO-MB2

γ -(6-Aminoethyl)-adenosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-833-MB2	40 μ l (1 mM)



Structural formula of γ -(6-Aminoethyl)-ATP-ATTO-MB2



absorption spectrum of ATTO MB2

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: C₃₅H₄₉N₉O₁₄P₃S (free acid)

Molecular Weight: 944.80 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{abs} 668 nm, ϵ 100.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)