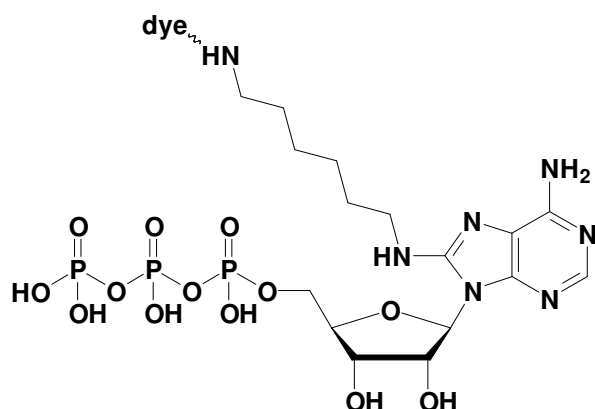


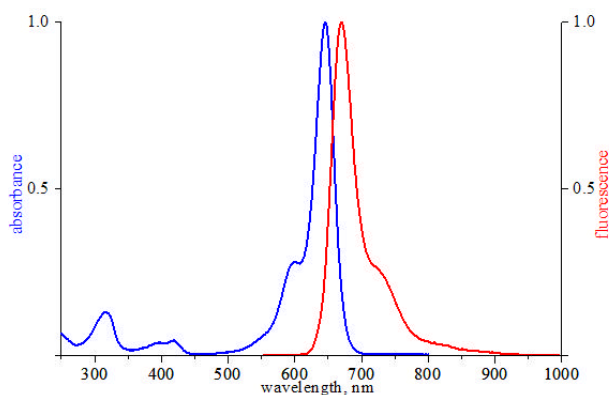
8-[(6-Amino)hexyl]-amino-ATP - ATTO 647N

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, labeled with ATTO 647N,
 Triethylammonium salt

Cat. No.	Amount
NU-807-647N	20 μ l / 1 mM



structural formula of 8-[(6-Amino)hexyl]-amino-ATP



excitation and emission spectrum of ATTO 647N

Cat. No.: NU-807-647N

Molecular Formula: C₁₆H₃₀N₇O₁₃P₃ - ATTO 647N
 (free acid)

Molecular Weight: 1248.37 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 644 nm; λ_{em} 669 nm; ϵ 150,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!