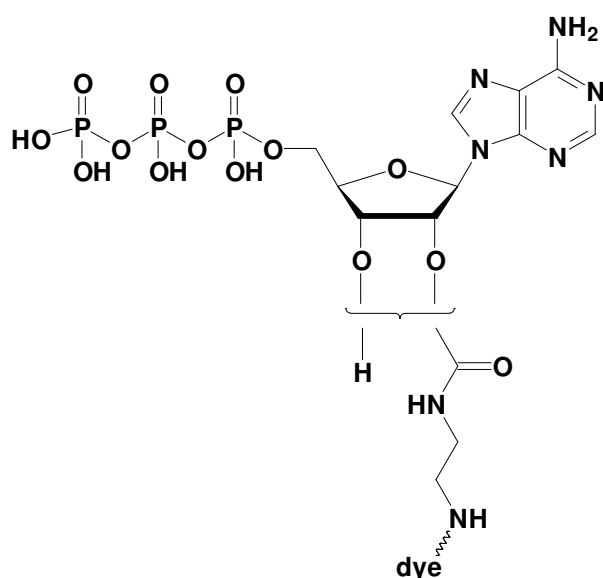


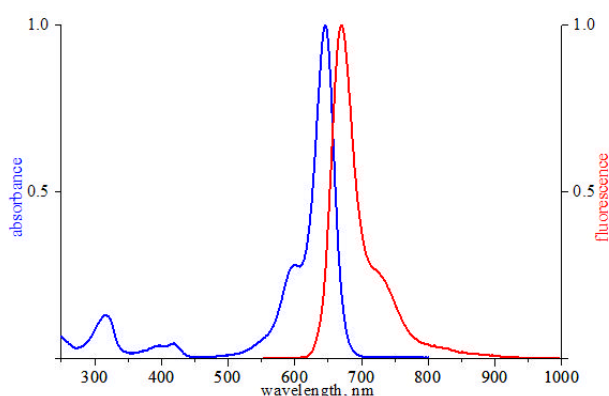
EDA-ATP - ATTO 647N

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-triphosphate, labeled with ATTO 647N,
 Triethylammonium salt

Cat. No.	Amount
NU-808-647N	10 μ l / 1 mM



structural formula of EDA-ATP



excitation and emission spectrum of ATTO 647N

Cat. No.: NU-808-647N

Molecular Formula: C₁₃H₂₂N₇O₁₄P₃ - ATTO 647N
 (free acid)

Molecular Weight: 1220.27 (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!

Selected References:

Hunke *et al.* (2010) The effect of NBD-Cl in nucleotide-binding of the major subunit alpha and B of the motor proteins F1FO ATP synthase and A1AO ATP synthase. *J. Bioenerg. Biomembr.* **42**:1.

Luo *et al.* (2008) Crystal structure of the NS3 protease-helicase from dengue virus. *J. Virol.* **82** (1):173.