



EDA-ADP-ATTO-647N

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

Cat. No.	Amount
NU-802-647N	40 µl (1 mM)

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: C₅₅H₇₀N₁₀O₁₃P₂ (free acid)

Molecular Weight: 1141.15 g/mol (free acid)

Exact Mass: 1140.46 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

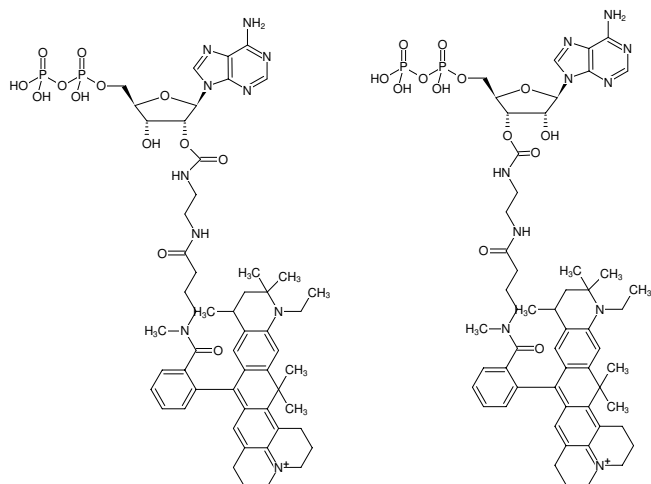
Form: solution in water

Color: blue

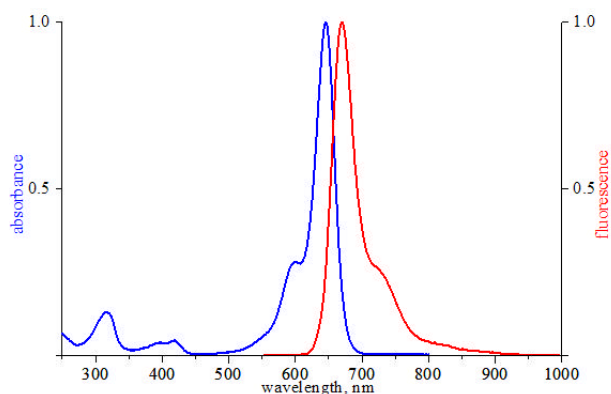
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ADP-ATTO-647N



excitation and emission spectrum of ATTO 647N

Selected References:

Singh *et al.* (2018) Crystallographic and enzymatic insights into the mechanisms of Mg-ADP inhibition in the A1 complex of the A1AO ATP synthase. *J. Struct. Biol.* **201** (1):26.

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):1.

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