



EDA-ADP-ATTO-495

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with ATTO 495, Triethylammonium salt

Cat. No.	Amount
NU-802-495	80 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: C₃₄H₄₄N₁₀O₁₂P₂ (free acid)

Molecular Weight: 846.72 g/mol (free acid)

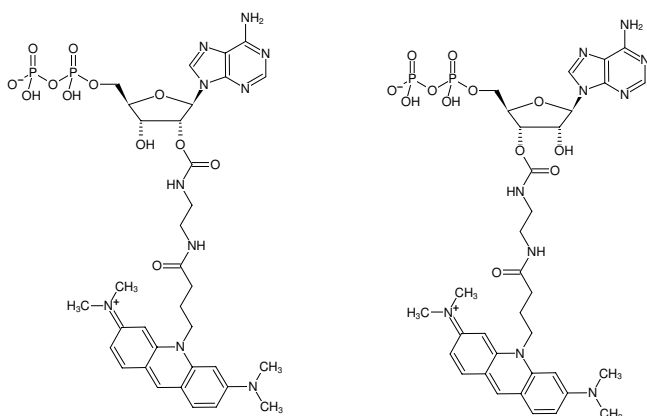
Purity: ≥ 95 % (HPLC)

Form: solution in water

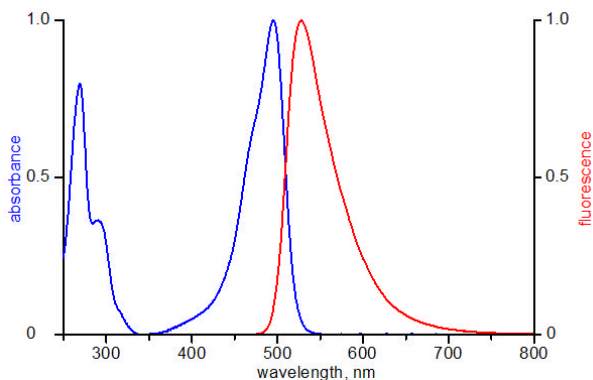
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 498 nm, λ_{em} 526 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ADP-ATTO-495



excitation and emission spectrum of ATTO 495

Selected References:

Wei *et al.* (2018) Adapted ATPase domain communication overcomes the cytotoxicity of p97 inhibitors. *J. Biol. Chem.* **293** (52):20169.