



EDA-ADP-ATTO-680

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with ATTO 680, Triethylammonium salt

Cat. No.	Amount
NU-802-680	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: C₄₀H₅₀N₁₀O₁₆P₂S (free acid)

Molecular Weight: 1020.90 g/mol (free acid)

Exact Mass: 1020.26 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

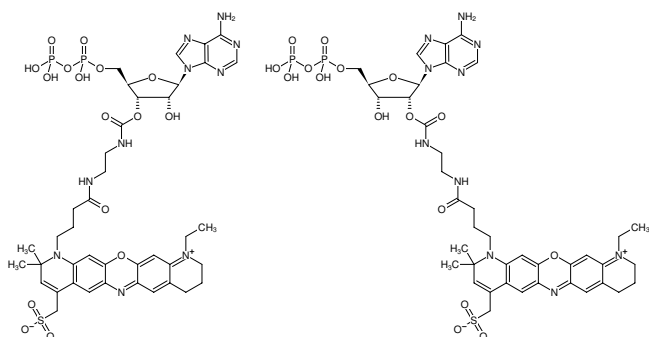
Form: solution in water

Color: blue

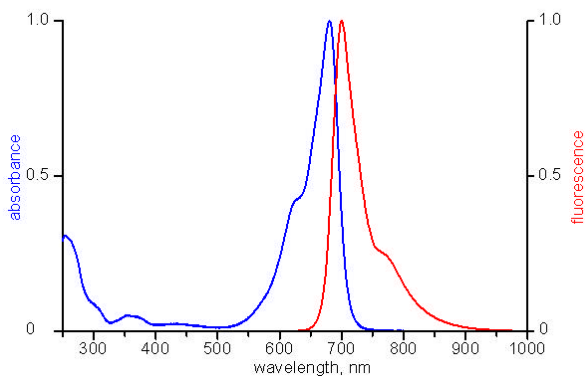
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 681 nm, λ_{em} 698 nm, ε 125.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-ADP-ATTO-680



excitation and emission spectrum of ATTO 680