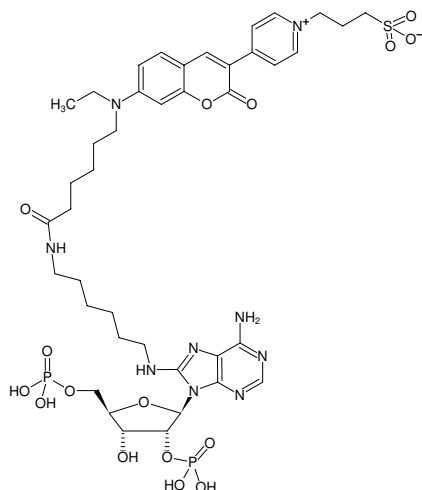




8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-DY-485XL

8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate, labeled with DY 485XL, Triethylammonium salt

Cat. No.	Amount
NU-812-485XL	40 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-adenosine-2',5'-bisphosphate-DY-485XL

For research use only!

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: C₄₁H₅₇N₉O₁₆P₂S (free acid)

Molecular Weight: 1025.96 g/mol (free acid)

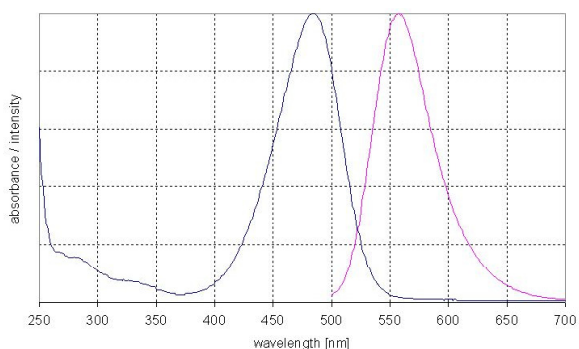
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 485 nm, λ_{em} 560 nm, ε 50.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of DY 485XL