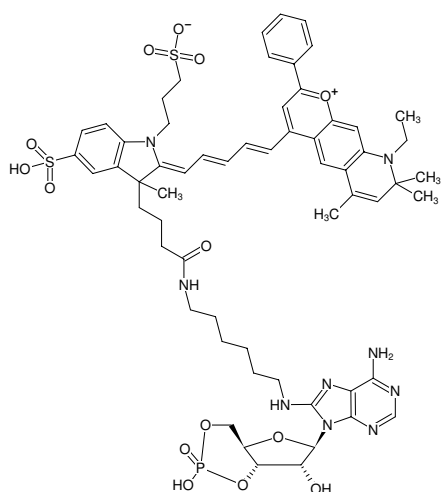




8-(6-Aminoethyl)-amino-cAMP-DY-776

8-(6-Aminoethyl)-amino-adenosine-3',5'-cyclic monophosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-851-776	80 µl (1 mM)



Structural formula of 8-(6-Aminoethyl)-amino-cAMP-DY-776

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: C₆₀H₇₂N₉O₁₄PS₂ (free acid)

Molecular Weight: 1238.37 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

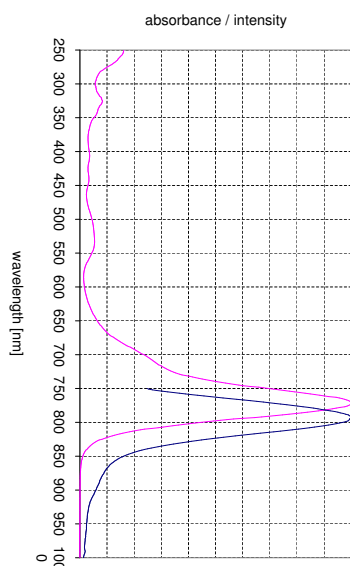
Form: solution in water

Color: dark blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 771 nm, λ_{em} 793 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of DY 776