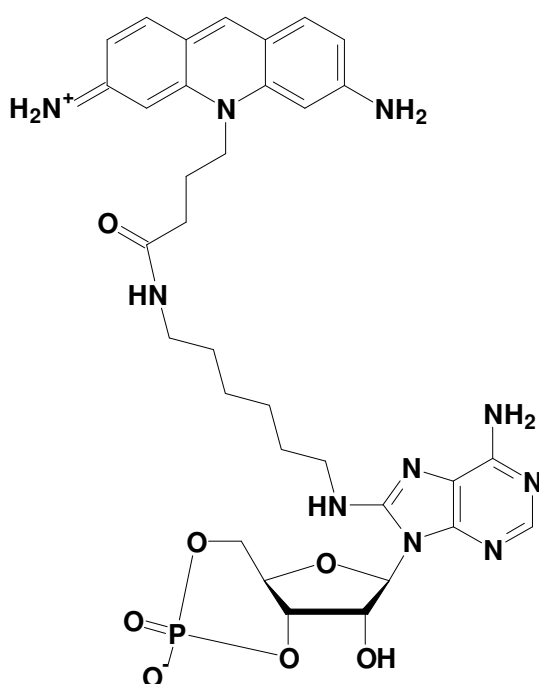


8-[(6-Amino)hexyl]-amino-cAMP - ATTO 465

8-[(6-Amino)hexyl]-amino-adenosine-3',5'-cyclic monophosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-851-465	120 μ l (1 mM)



structural formula of 8-[(6-Amino)hexyl]-amino-cAMP - ATTO 465

Cat. No.: NU-851-465

Molecular Formula: C₃₃H₄₁N₁₀O₇P (free acid)

Molecular Weight: 720.72 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

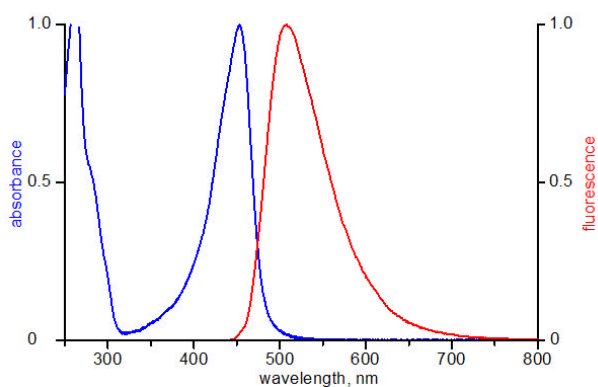
Spectroscopic properties:

λ_{exc} 453 nm; λ_{em} 508 nm; ϵ 75,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!



excitation and emission spectrum of ATTO 465