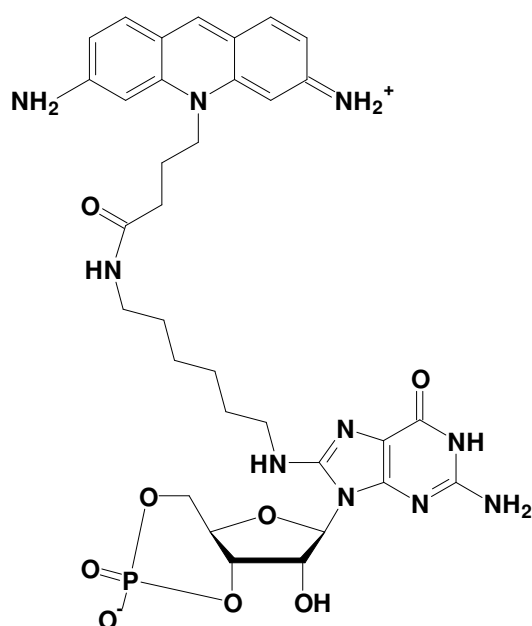


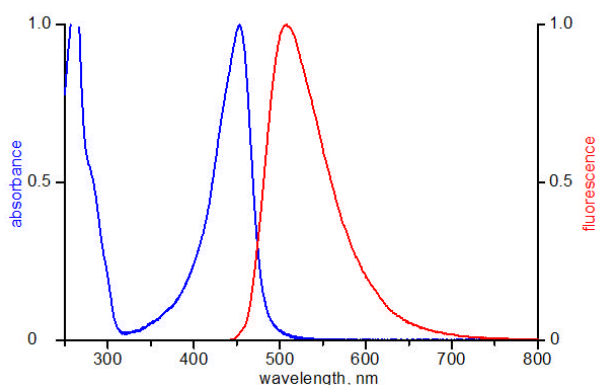
8-[(6-Amino)hexyl]-amino-cGMP - ATTO 465

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

Cat. No.	Amount
NU-832-465	120 μ l (1 mM)



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



excitation and emission spectrum of ATTO 465

Cat. No.: NU-832-465

Molecular Formula: C₃₃H₄₁N₁₀O₈P (free acid)

Molecular Weight: 736.72 (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!