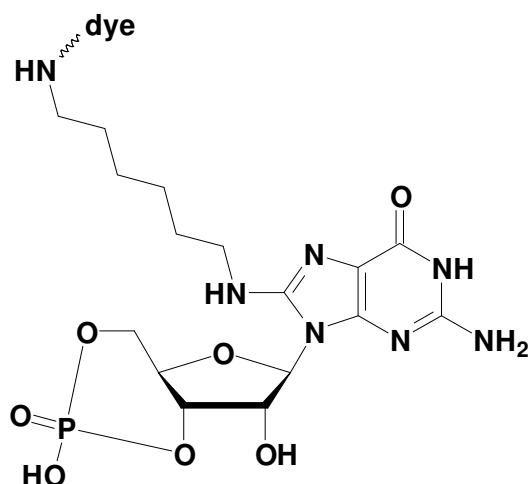


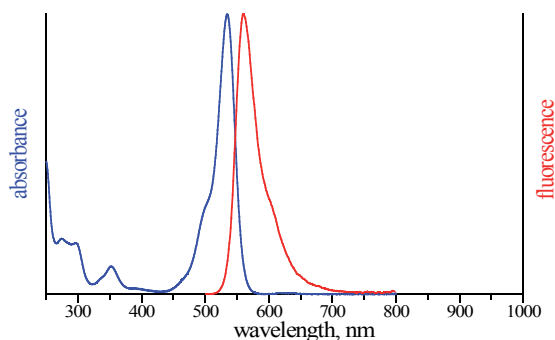
8-[(6-Amino)hexyl]-amino-cGMP - ATTO Rho6G

8-[(6-Amino)hexyl]-amino-guanosine-3',5'-cyclic monophosphate, labeled with ATTO Rho6G, Triethylammonium salt

Cat. No.	Amount
NU-832-RHO6G	80 μ l (1 mM)



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



excitation and emission spectrum of ATTO Rho6G

Cat. No.: NU-832-RHO6G

Molecular Formula: C₁₆H₂₆N₇O₇P - ATTO Rho6G (free acid)

Molecular Weight: 954,39 (free acid)

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

Spectroscopic properties:

λ_{exc} 535 nm; λ_{em} 560 nm; ϵ 1,15 x 10⁵ 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!