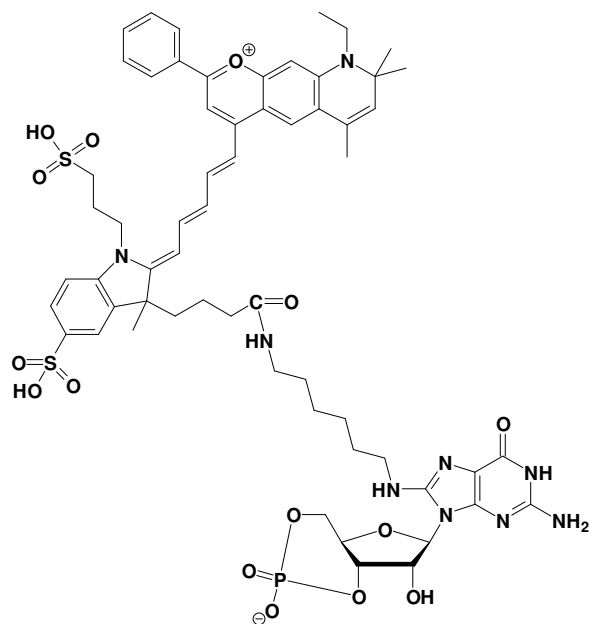


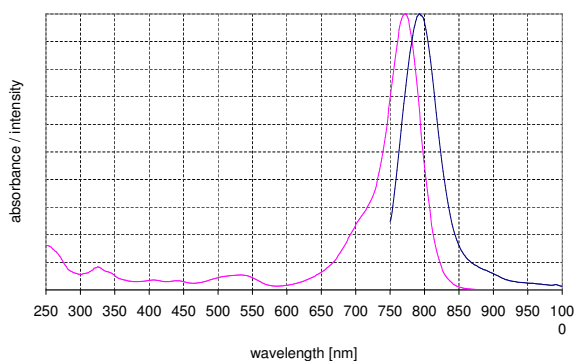
8-[(6-Amino)hexyl]-amino-cGMP - DY 776

8-[(6-Amino)hexyl]-amino-guanosine-3',5'-cyclic monophosphate, labeled with DY 776

Cat. No.	Amount
NU-832-776	20 μ l / 1 mM



structural formula of 8-[(6-Amino)hexyl]-amino-cGMP - DY 776



excitation and emission spectrum of DY 776

Cat. No.: NU-832-776

Molecular Formula: C₆₀H₇₂N₉O₁₅PS₂ (Anion)

Molecular Weight: 1254.39 (Anion)

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

Spectroscopic properties:

λ_{exc} 771 nm; λ_{em} 801 nm; ϵ 120,000 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!