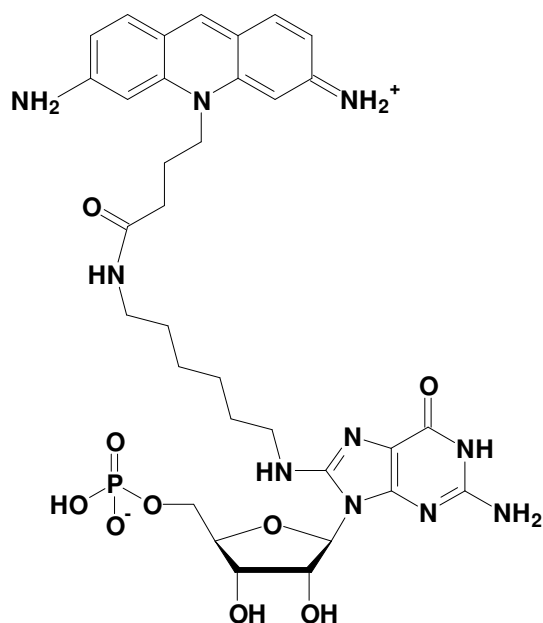


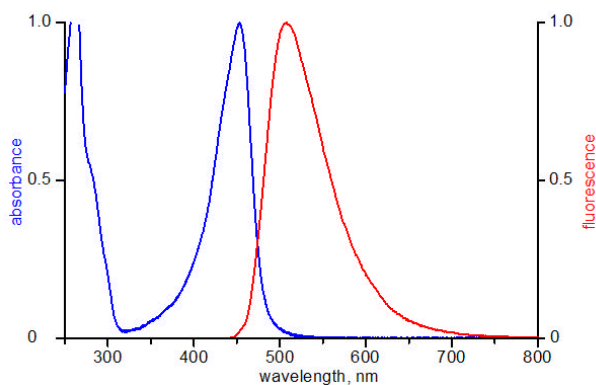
8-[(6-Amino)hexyl]-amino-GMP - ATTO 465

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

| Cat. No. | Amount |
|------------|--------------------|
| NU-829-465 | 120 μ l (1 mM) |



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



excitation and emission spectrum of ATTO 465

Cat. No.: NU-829-465

Molecular Formula: C₃₃H₄₃N₁₀O₉P (free acid)

Molecular Weight: 754.73 (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!