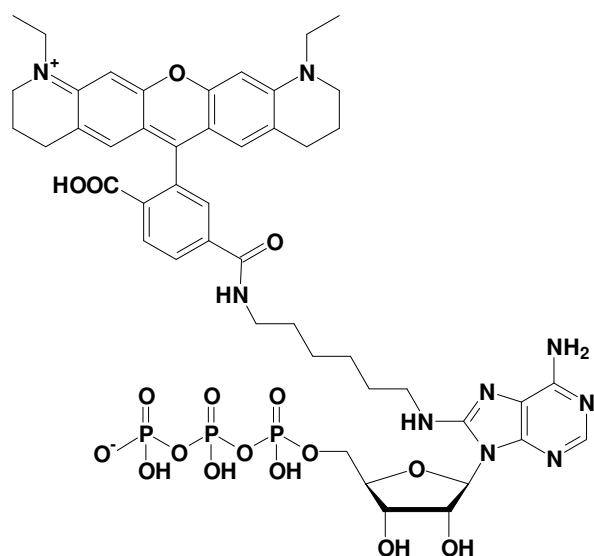


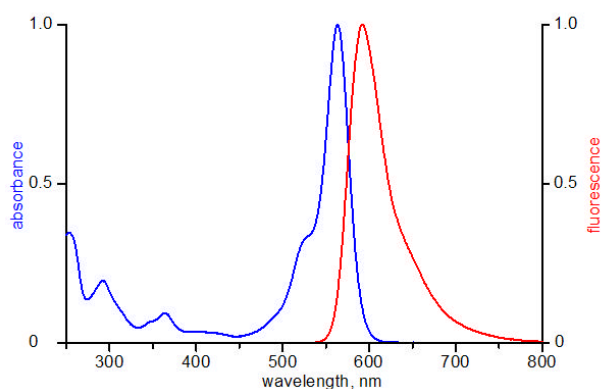
8-[(6-Amino)hexyl]-amino-ATP - ATTO 565

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, labeled with ATTO 565,
 Triethylammonium salt

Cat. No.	Amount
NU-807-565	30 μ l / 1 mM



structural formula of 8-[(6-Amino)hexyl]-amino-ATP - ATTO 565



excitation and emission spectrum of ATTO 565

Cat. No.: NU-807-565

Molecular Formula: C₄₇H₅₈N₉O₁₇P₃ (free acid)

Molecular Weight: 1113.94 (free acid)

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

Spectroscopic properties:

λ_{exc} 563 nm; λ_{em} 592 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!