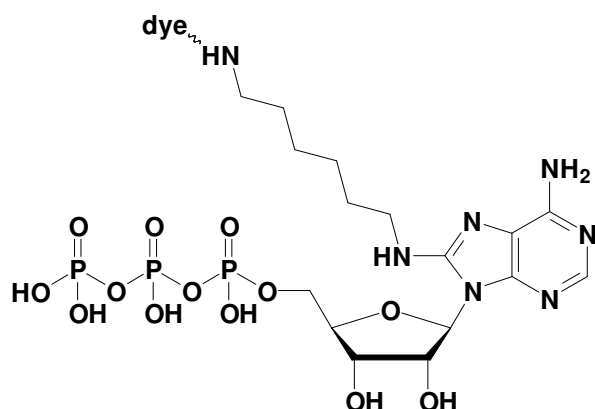


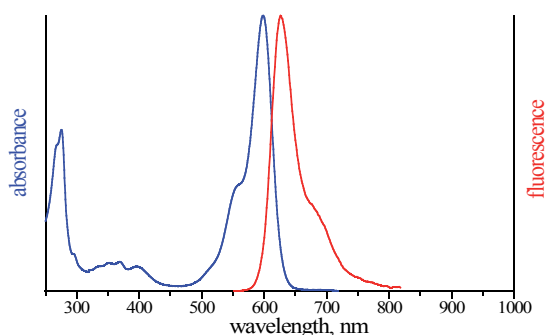
8-[(6-Amino)hexyl]-amino-ATP - ATTO Rho13

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, labeled with ATTO Rho13,
 Triethylammonium salt

Cat. No.	Amount
NU-807-RHO13	80 μ l (1 mM)



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



excitation and emission spectrum of ATTO Rho13

Cat. No.: NU-807-RHO13

Molecular Formula: C₁₆H₃₀N₇O₁₃P₃ - ATTO Rho13
 (free acid)

Molecular Weight: 1248,37 (free acid)

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

Spectroscopic properties:

λ_{exc} 600 nm; λ_{em} 625 nm; ϵ 1,25 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!