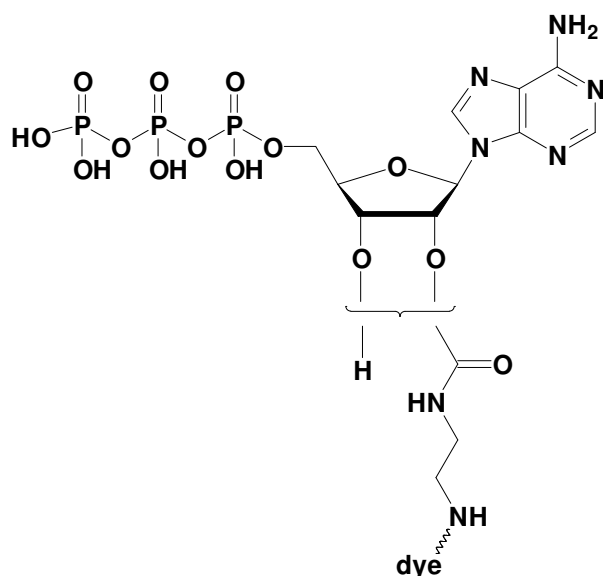


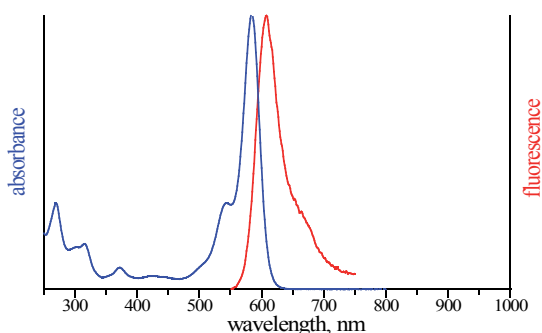
EDA-ATP - ATTO Rho101

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, labeled with ATTO Rho101, Triethylammonium salt

Cat. No.	Amount
NU-808-RHO101	50 μ l / (0.2 mM)



structural formula of EDA-ATP



excitation and emission spectrum of ATTO Rho101

Cat. No.: NU-808-RHO101

Molecular Formula: C₁₃H₂₂N₇O₁₄P₃ - ATTO Rho101 (free acid)

Molecular Weight: 1164,27 (free acid)

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

Spectroscopic properties:

λ_{exc} 586 nm; λ_{em} 610 nm; ϵ 1,2 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!