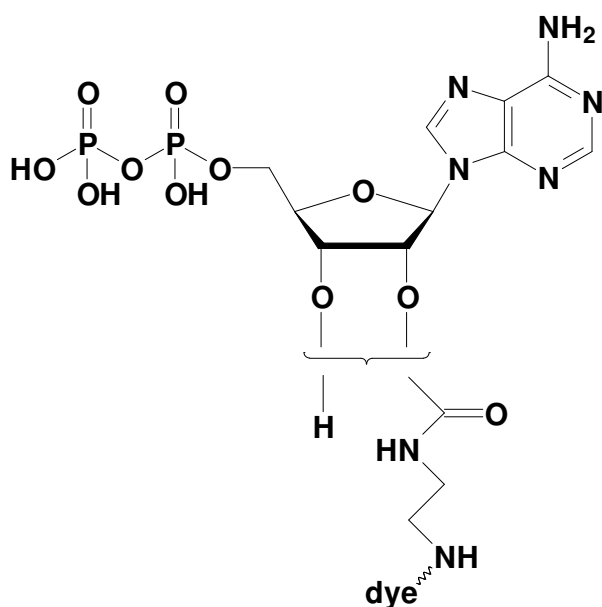


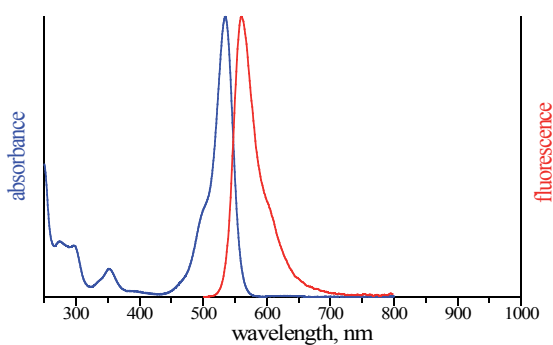
EDA-ADP - ATTO Rho6G

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with ATTO Rho6G, Triethylammonium salt

Cat. No.	Amount
NU-802-RHO6G	40 μ l (1 mM)



structural formula of EDA-ADP



excitation and emission spectrum of ATTO Rho6G

Cat. No.: NU-802-RHO6G

Molecular Formula: C₁₃H₂₁N₇O₁₁P₂ - ATTO Rho6G (free acid)

Molecular Weight: 1008,29 (Anion)

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

Spectroscopic properties:

λ_{exc} 535 nm; λ_{em} 560 nm; ϵ 1,15 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!