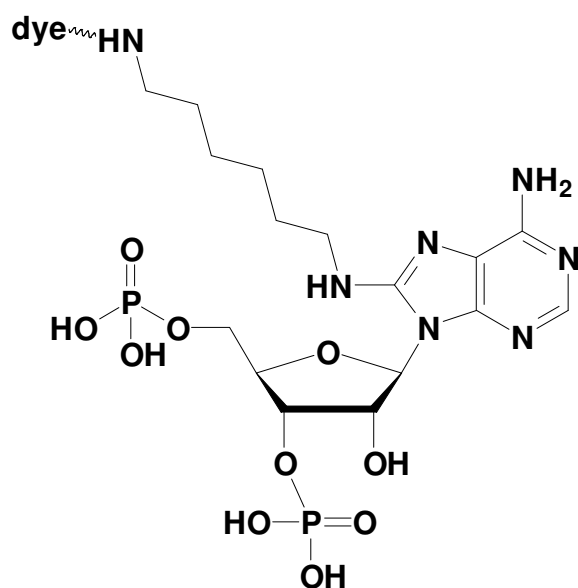


8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate - ATTO Rho6G 8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate, labeled with ATTO Rho6G, Triethylammonium salt

Cat. No.	Amount
NU-811-RHO6G	40 μ l (1 mM)



Cat. No.: NU-811-RHO6G

Molecular Formula: C₁₆H₂₉N₇O₁₀P₃ - ATTO Rho6G
 (free acid)

Molecular Weight: 1036,39 (free acid)

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!

