

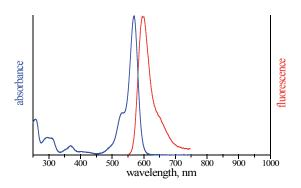


8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho11

8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-812-RHO11	40 μl (1 mM)

Structural formula of 8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient

temperature possible.

Shelf Life: 12 months after date of delivery **Molecular Formula:** C₅₁H₆₆N₁₀O₁₃P₂ (free acid) **Molecular Weight:** 1089.09 g/mol (free acid)

Exact Mass: 1088.43 g/mol (free acid)

Purity: ≥ 95 % (HPLC) **Form:** solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 572 nm, λ_{em} 595 nm,

 ϵ 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)