

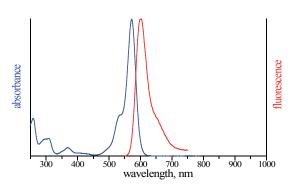


8-(6-Aminohexyl)-amino-adenosine-3',5'-bisphosphate-ATTO-Rho12

8-(6-Aminohexyl)-amino-adenosine-3',5'-bisphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-811-RHO12	40 μl (1 mM)

Structural formula of 8-(6-Aminohexyl)-amino-adenosine-3',5'-bisphosphate-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12

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: $C_{57}H_{78}N_{10}O_{13}P_2$ (free acid) Molecular Weight: 1173.26 g/mol (free acid)

Exact Mass: 1172.52 g/mol (free acid)

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

Concentration: 1.0 mM - 1.1 mM

Spectroscopic Properties: λ_{exc} 576 nm, λ_{em} 601 nm,

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