

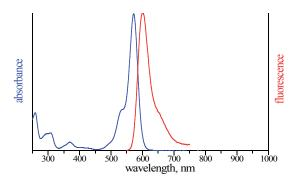


y-(6-Aminohexyl)-ATP-ATTO-Rho12

γ-(6-Aminohexyl)-adenosine-5'-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-833-RHO12	80 μl (1 mM)

Structural formula of γ -(6-Aminohexyl)-ATP-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:** $C_{57}H_{78}N_9O_{16}P_3$ (free acid) **Molecular Weight:** 1238.22 g/mol (free acid)

Exact Mass: 1237.48 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} 576 nm, λ_{em} 601 nm,

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