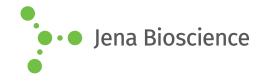
DATA SHEET



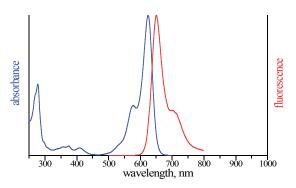


EDA-ATP-ATTO-Rho14

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, labeled with ATTO Rho14, Triethylammonium salt

Cat. No.	Amount
NU-808-RHO14	40 μl (1 mM)

Structural formula of EDA-ATP-ATTO-Rho14



excitation and emission spectrum of ATTO Rho14

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:** $C_{54}H_{63}Cl_4N_{10}O_{17}P_3$ (free acid) **Molecular Weight:** 1358.87 g/mol (free acid)

Exact Mass: 1356.23 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} 626 nm, λ_{em} 646 nm,

ε 140.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)