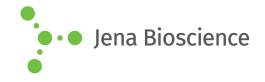
DATA SHEET

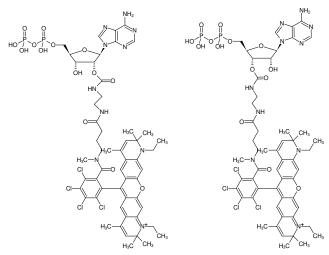




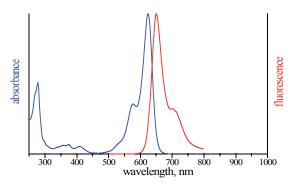
EDA-ADP-ATTO-Rho14

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

Cat. No.	Amount
NU-802-RHO14	40 μl (1 mM)



Structural formula of EDA-ADP-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: $C_{54}H_{62}Cl_4N_{10}O_{14}P_2$ (free acid) Molecular Weight: 1278.89 g/mol (free acid)

Exact Mass: 1276.27 g/mol (free acid)

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

Color: blue

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)