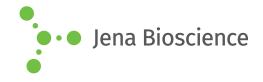
DATA SHEET

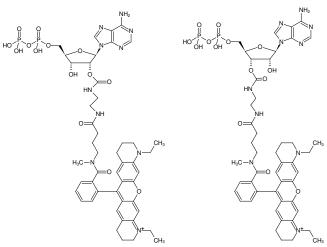




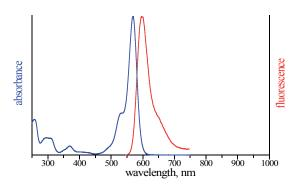
EDA-ADP-ATTO-Rho11

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-802-RHO11	40 μl (1 mM)



Structural formula of EDA-ADP-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:** $C_{48}H_{58}N_{10}O_{14}P_2$ (free acid) **Molecular Weight:** 1061.00 g/mol (free acid)

Exact Mass: 1060.36 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)