

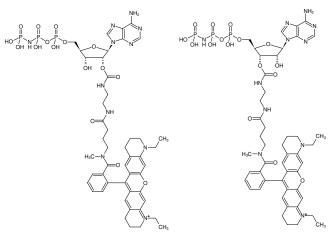




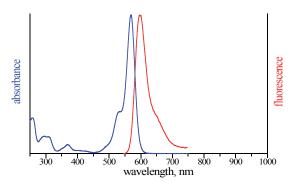
## EDA-AppNHp (EDA-AMPPNP)-ATTO-Rho11

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-[(β,γ)-imido] triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-810-RHO11	40 µl (1 mM)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-ATTO-Rho11



excitation and emission spectrum of ATTO Rho11

## For general laboratory use.

Shipping: shipped on dry ice Storage Conditions: store at -20 °C Shelf Life: 6 months after date of delivery Molecular Formula:  $C_{48}H_{60}N_{11}O_{16}P_3$  (free acid) Molecular Weight: 1139.99 g/mol (free acid) Exact Mass: 1139.34 g/mol (free acid) Purity:  $\ge 95$  % (HPLC) Form: solution in water Concentration: 1.0 mM - 1.1 mM

## **pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  572 nm,  $\lambda_{em}$  595 nm,  $\epsilon$  120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



rtified QMS and EMS according to N EN ISO 9001 and DIN EN ISO 14001 g.-No.: ICV03597 034 and ICV03597 534