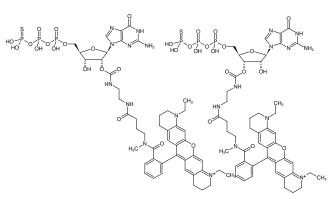




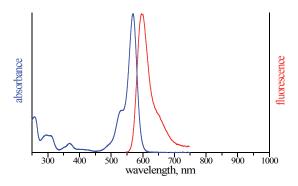
## EDA-GTPyS-ATTO-Rho11

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO Rho11, Triethylammonium salt

Cat. No.	Amount
NU-1610-RHO11	50 µl (1 mM)



Structural formula of EDA-GTPyS-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:  $C_{48}H_{59}N_{10}O_{17}P_3S$  (free acid) Molecular Weight: 1173.03 g/mol (free acid) Exact Mass: 1172.30 g/mol (free acid) Purity:  $\ge 90$  % (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)

