



### EDA-GTPyS-ATTO-Rho12

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate, labeled with ATTO Rho12, Triethylammonium salt

Cat. No.	Amount
NU-1610-RHO12	50 µl (1 mM)

**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<sub>54</sub>H<sub>71</sub>N<sub>10</sub>O<sub>17</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1257.20 g/mol (free acid)

**Exact Mass:** 1256.39 g/mol (free acid)

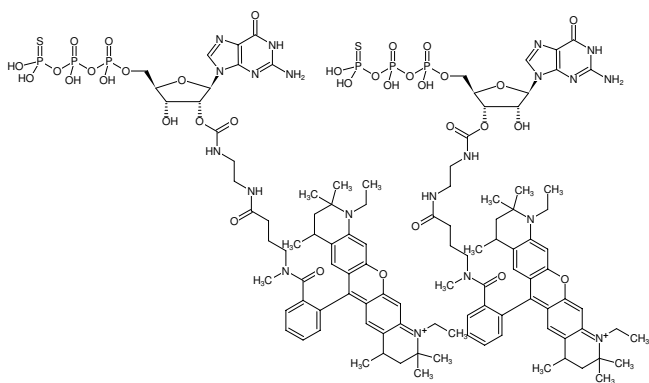
**Purity:** ≥ 90 % (HPLC)

**Form:** solution in water

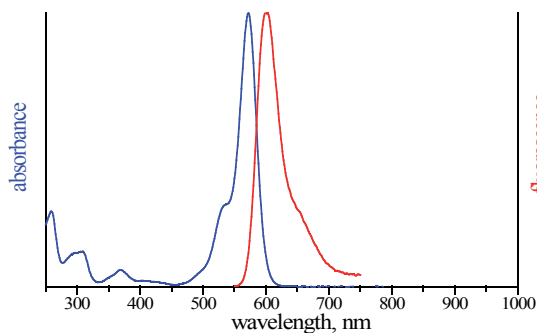
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 576 nm, λ<sub>em</sub> 601 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-GTPyS-ATTO-Rho12



excitation and emission spectrum of ATTO Rho12