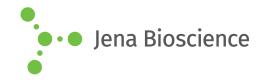
DATA SHEET



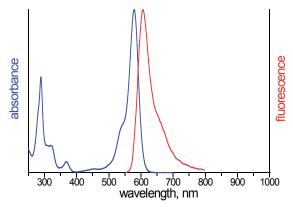


EDA-GTPyS-ATTO-Thio12

 $2^\prime/3^\prime\text{-O-}(2\text{-Aminoethyl-carbamoyl})\text{-guanosine-5}^\prime\text{-}(\gamma\text{-thio})\text{-triphosphate, labeled with ATTO Thio12, Triethylammonium salt}$

Cat. No.	Amount
NU-1610-THIO12	50 μl (1 mM)

Structural formula of EDA-GTPyS-ATTO-Thio12



excitation and emission spectrum of ATTO Thio12

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: $C_{42}H_{51}N_{10}O_{16}P_3S_2$ (free acid)

Molecular Weight: 1108.96 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_{exc} 582 nm, λ_{em} 607 nm, ϵ 110.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)