



EDA-GDP-ATTO-532

2'/3'-O-(2-Aminoethyl-carbamoyl)-Guanosine-5'-diphosphate, labeled with ATTO 532, Triethylammonium salt

Cat. No.	Amount
NU-840-532	80 µl (1 mM)

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: C₄₂H₅₀N₁₀O₂₁P₂S₂ (free acid)

Molecular Weight: 1156.98 g/mol (free acid)

Exact Mass: 1156.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

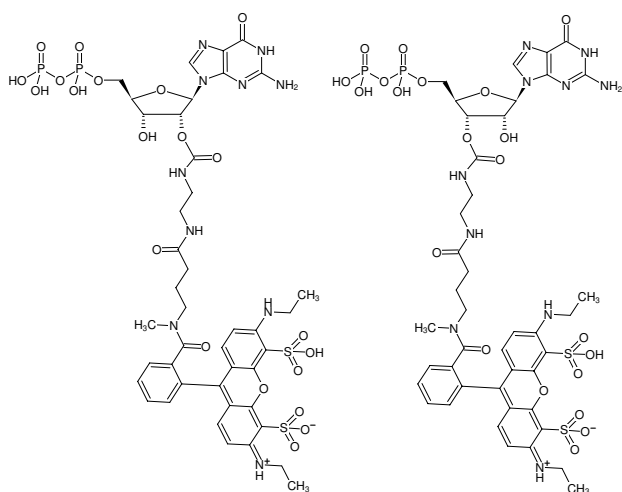
Form: solution in water

Color: orange

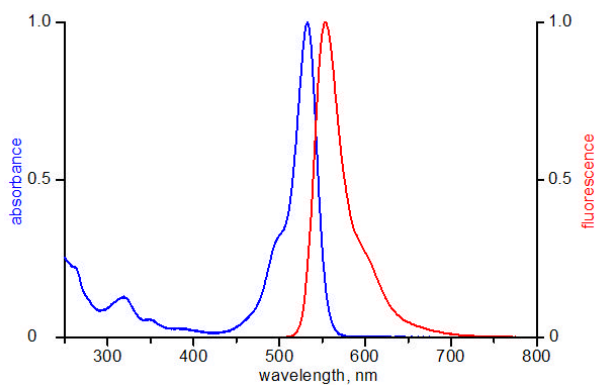
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 532 nm, λ_{em} 552 nm, ε 115.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-GDP-ATTO-532



excitation and emission spectrum of ATTO 532