

**EDA-m⁷GDP-ATTO-550**

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 550, Triethylammonium salt

Cat. No.	Amount
NU-827-550	80 µl (1 mM)

For research use only!

Shipping: shipped on blue ice

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: C₅₁H₆₄N₁₀O₁₅P₂ (free acid)

Molecular Weight: 1119.08 g/mol (free acid)

Exact Mass: 1118.40 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

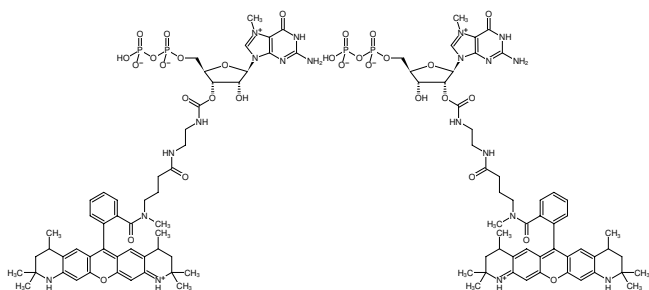
Form: solution in water

Color: red-violet

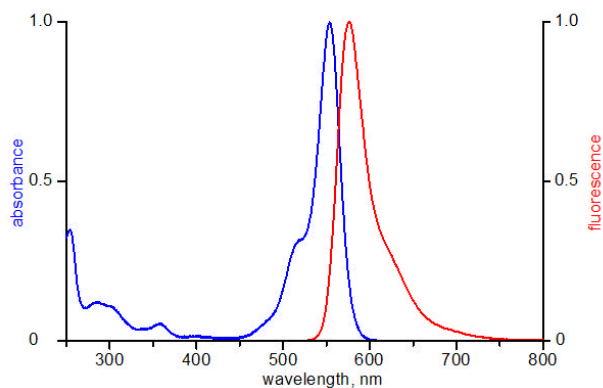
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 554 nm, λ_{em} 576 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-ATTO-550



excitation and emission spectrum of ATTO 550