



EDA-m⁷GDP-ATTO-590

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 590

Cat. No.	Amount
NU-827-590	120 µl (1 mM)

For research use only!

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: C₅₁H₅₉N₉O₁₆P₂ (free acid)

Molecular Weight: 1116.02 g/mol (free acid)

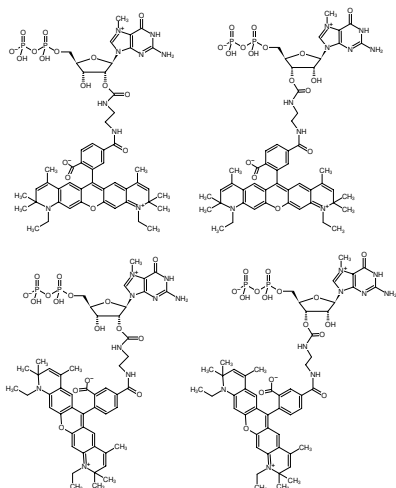
Purity: ≥ 95 % (HPLC)

Form: solution in water

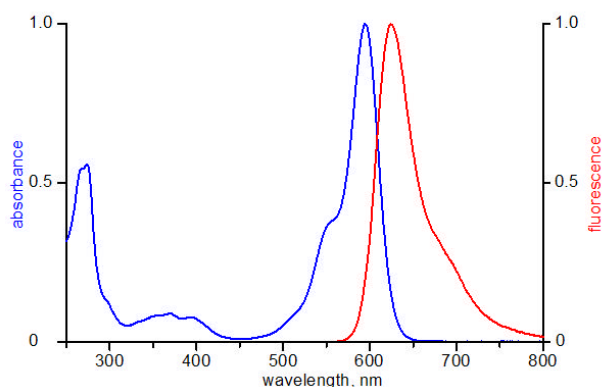
Concentration: 1.0 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 593 nm, λ_{em} 622 nm, ε 120.0 L mol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-ATTO-590



excitation and emission spectrum of ATTO 590