



EDA-m⁷GDP-6-JOE

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with 6-JOE, Triethylammonium salt

Cat. No.	Amount
NU-827-JOE	160 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: C₃₇H₃₅Cl₂N₇O₂₀P₂ (free acid)

Molecular Weight: 1029.08 g/mol (free acid)

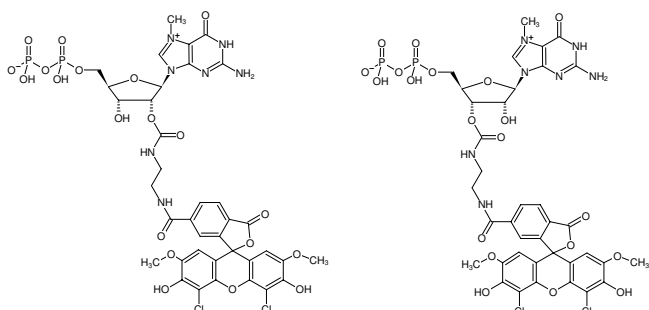
Purity: ≥ 95 % (HPLC)

Form: solution in water

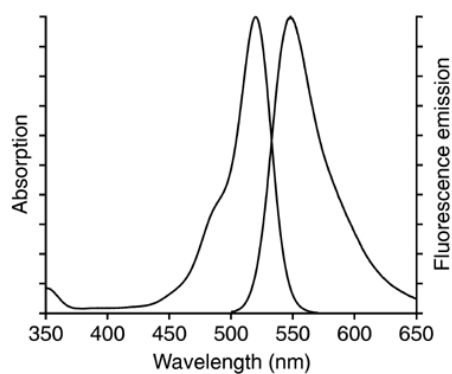
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 520 nm, λ_{em} 548 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of EDA-m⁷GDP-6-JOE



excitation and emission spectrum of 6-JOE