



7-Propargylamino-7-deaza-dGTP-6-JOE

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with 6-JOE, Triethylammonium salt

Cat. No.	Amount
NU-1615-JOE	120 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: C₃₇H₃₂Cl₂N₅O₂₁P₃ (free acid)

Molecular Weight: 1045.02 g/mol (free acid)

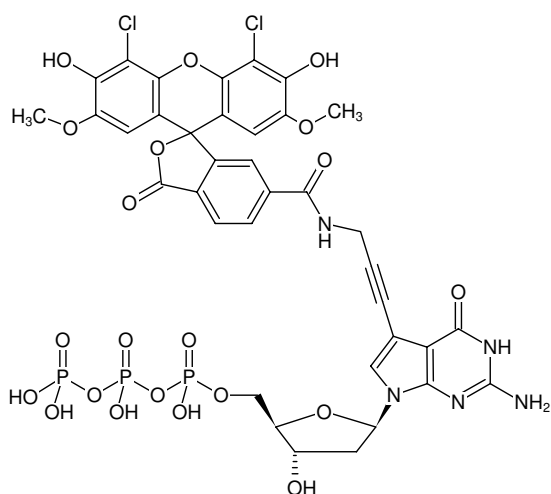
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

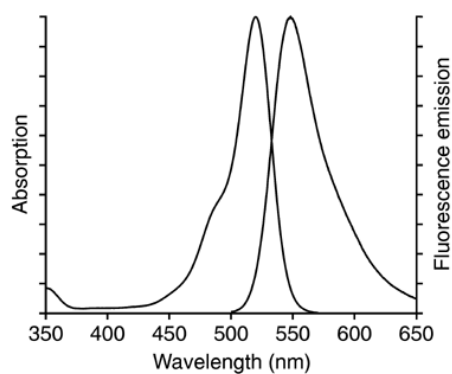
Concentration: 1.0 mM - 1.1 mM

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 7-Propargylamino-7-deaza-dGTP-6-JOE



excitation and emission spectrum of 6-JOE