



7-Propargylamino-7-deaza-ddGTP-DY-751

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with DY 751, Triethylammonium salt

| Cat. No. | Amount |
|-------------|----------------|
| NU-1618-751 | 80 µl (0,5 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: C₅₆H₇₀N₇O₂₀P₃S₂ (free acid)

Molecular Weight: 1317.33 g/mol (free acid)

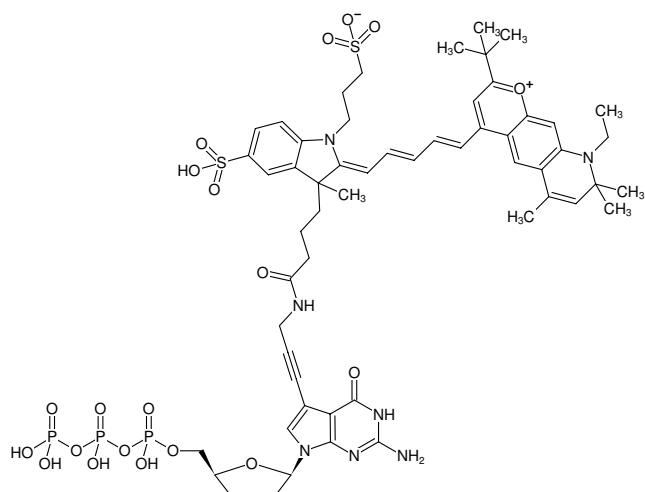
Purity: ≥ 95 % (HPLC)

Form: solution in water

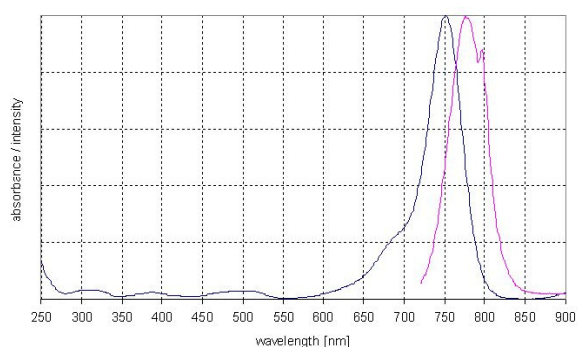
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 751 nm, λ_{em} 779 nm, ε 270.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of 7-Propargylamino-7-deaza-ddGTP-DY-751



excitation and emission spectrum of DY 751