



5-Propargylamino-ddCTP-DY-751

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with DY 751, Triethylammonium salt

| Cat. No. | Amount |
|------------|----------------|
| NU-850-751 | 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: C₅₅H₇₀N₅O₂₀P₃S₂ (free acid)

Molecular Weight: 1278.22 g/mol (free acid)

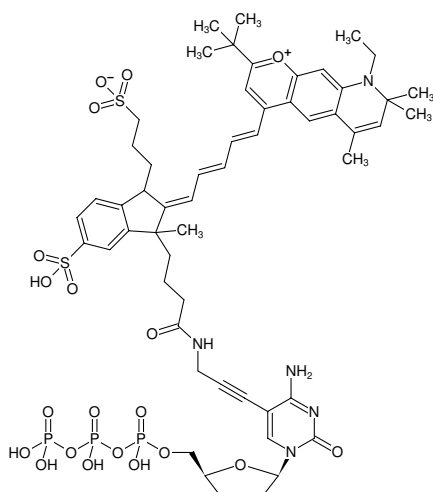
Purity: ≥ 95 % (HPLC)

Form: solution in water

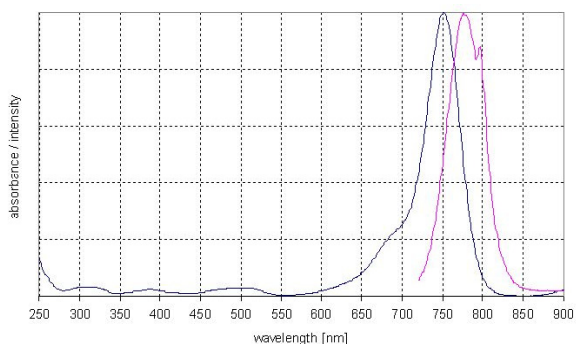
Concentration: 0.50 mM - 0.55 mM

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 5-Propargylamino-ddCTP-DY-751



excitation and emission spectrum of DY 751