



5-Propargylamino-dCTP-ATTO-665

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with ATTO 665, Triethylammonium salt

Cat. No.	Amount
NU-809-665	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: C₁₂H₁₆N₄O₁₃P₃ - ATTO 665 (free acid)

Molecular Weight: 1124.22 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

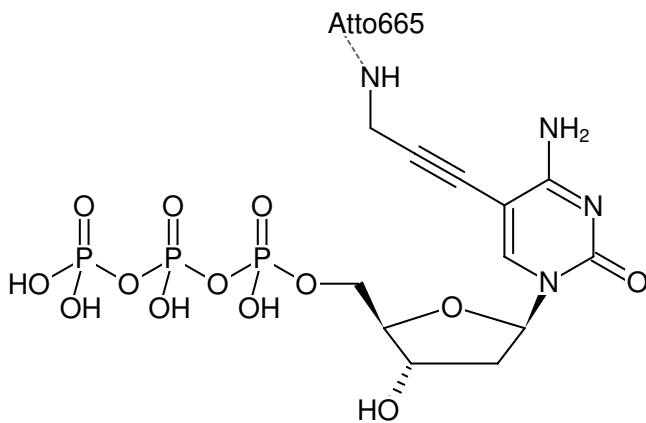
Form: solution in water

Color: blue

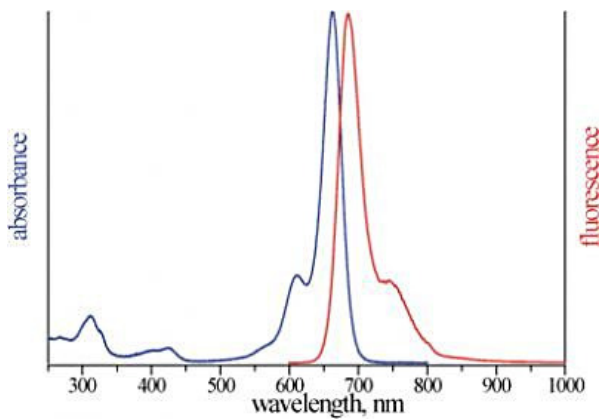
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 662 nm, λ_{em} 680 nm, ε 160.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 5-Propargylamino-dCTP-ATTO-665



excitation and emission spectrum of ATTO 665