



7-Propargylamino-7-deaza-ddGTP-ATTO-665

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with ATTO 665, Triethylammonium salt

Cat. No.	Amount
NU-1618-665	400 µl (0,1 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: C₁₄H₂₀N₅O₁₂P₃ - ATTO 665 (free acid)

Molecular Weight: 1147.26 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

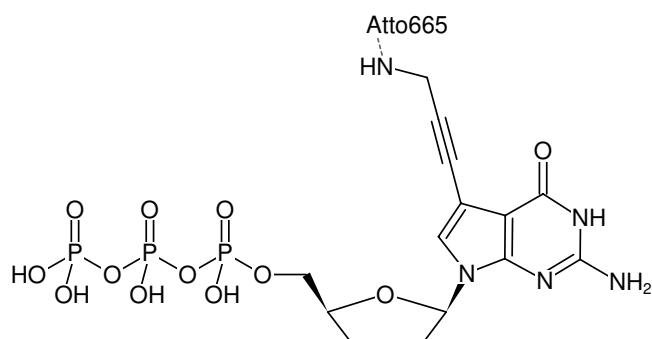
Form: solution in water

Color: blue

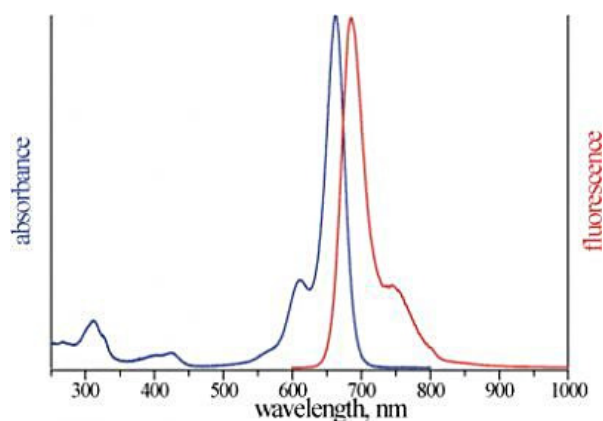
Concentration: 0.10 mM - 0.11 mM

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 7-Propargylamino-7-deaza-ddGTP-ATTO-665



excitation and emission spectrum of ATTO 665