



7-Propargylamino-7-deaza-ddGTP-ATTO-740

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with ATTO 740, Triethylammonium salt

Cat. No.	Amount
NU-1618-740	80 µl (0,5 mM)

For research use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: long time storage at pH 7.5

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₄H₂₀N₅O₁₂P₃ - ATTO 740 (free acid)

Molecular Weight: 992.26 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

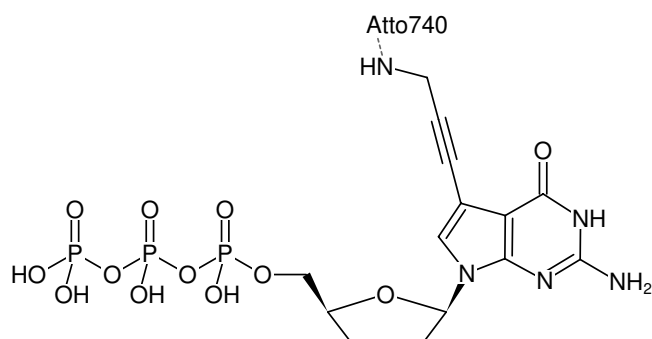
Form: solution in water

Concentration: 0.50 mM - 0.55 mM

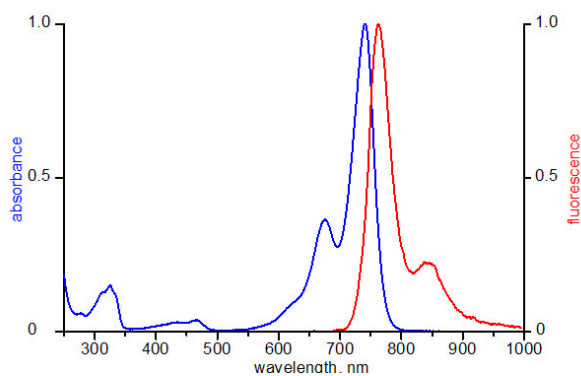
pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 743 nm, λ_{em} 763 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Please note: The dye decomposes at pH-levels greater 8.0. The recommended pH for a solution of this product is 7.5.



Structural formula of 7-Propargylamino-7-deaza-ddGTP-ATTO-740



excitation and emission spectrum of ATTO 740