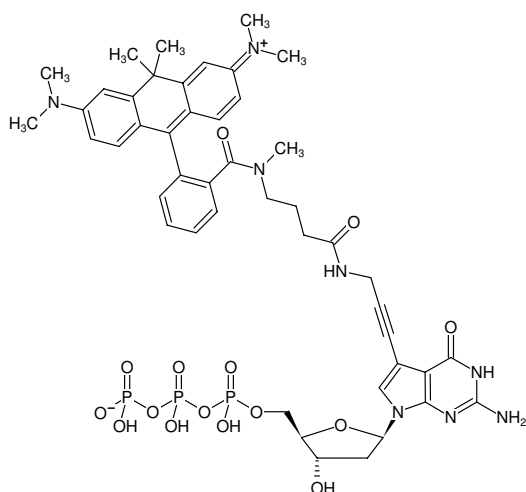




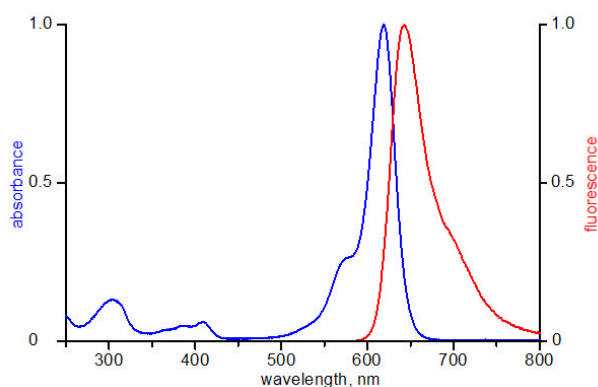
7-Propargylamino-7-deaza-dGTP-ATTO-620

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with ATTO 620, Triethylammonium salt

Cat. No.	Amount
NU-1615-620	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dGTP-ATTO-620



excitation and emission spectrum of ATTO 620

For research use only!

Shipping: shipped on blue ice

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: C₄₆H₅₅N₈O₁₅P₃ (free acid)

Molecular Weight: 1052.91 g/mol (free acid)

Exact Mass: 1052.30 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 620 nm, λ_{em} 642 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)