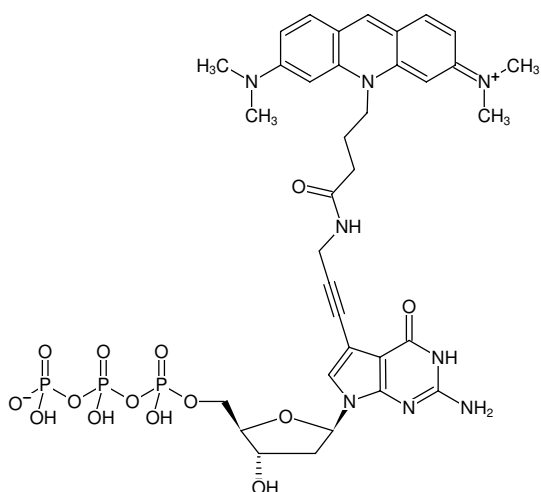




7-Propargylamino-7-deaza-ATTO-495

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate-, labeled with ATTO 495, Triethylammonium salt

Cat. No.	Amount
NU-1615-495	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dGTP-ATTO-495

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: C₃₅H₄₃N₈O₁₄P₃ (free acid)

Molecular Weight: 892.68 g/mol (free acid)

Exact Mass: 892.21 g/mol (free acid)

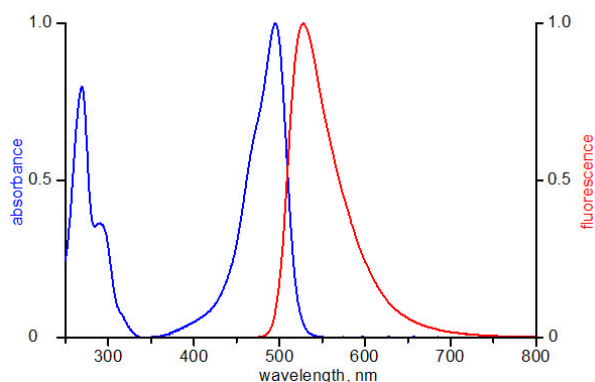
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 498 nm, λ_{em} 526 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 495