



7-Propargylamino-7-deaza-ddGTP-ATTO-425

7-Deaza-7-propargylamino-2',3'-dideoxyguanosine-5'-triphosphate, labeled with ATTO 425, Triethylammonium salt

Cat. No.	Amount
NU-1618-425	160 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: C₃₆H₄₅N₆O₁₇P₃ (free acid)

Molecular Weight: 926.70 g/mol (free acid)

Exact Mass: 926.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

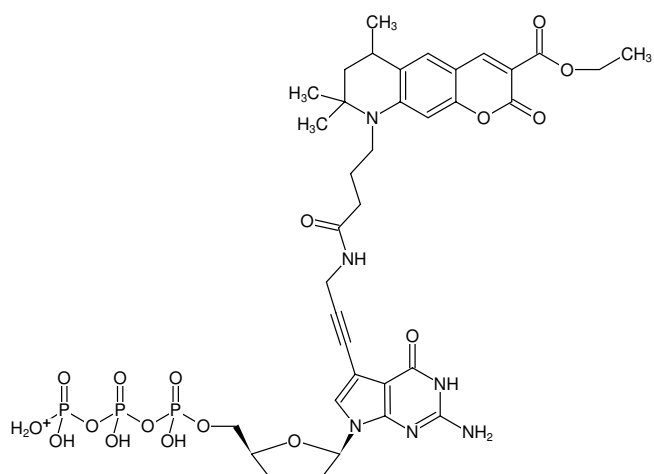
Form: solution in water

Color: green-yellow

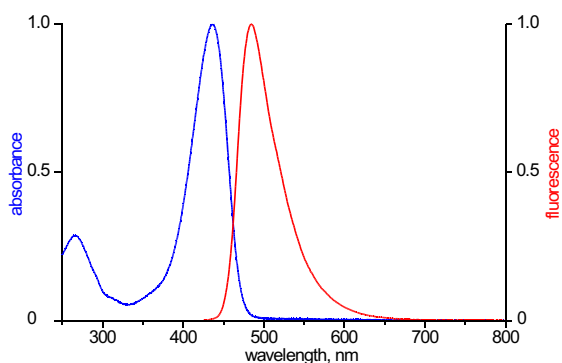
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 439 nm, λ_{em} 485 nm, ε 45.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ddGTP-ATTO-425



excitation and emission spectrum of ATTO 425