



7-Propargylamino-7-deaza-ddATP-ATTO-465

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with ATTO 465, Triethylammonium salt

Cat. No.	Amount
NU-1612-465	160 µl (0,5 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 2 ml).

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: C₃₁H₃₅N₈O₁₂P₃ (free acid)

Molecular Weight: 804.58 g/mol (free acid)

Exact Mass: 804.16 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

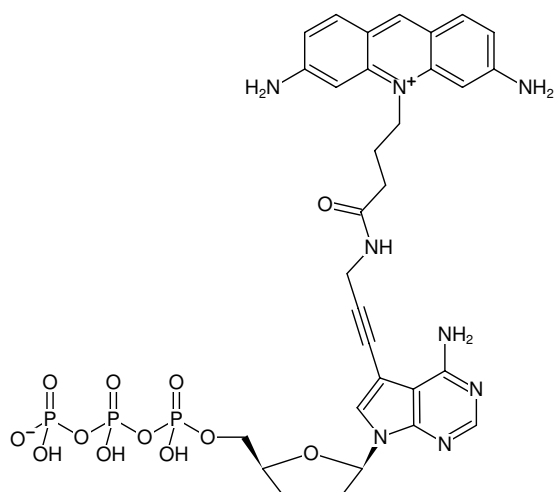
Form: solution in water

Color: yellow

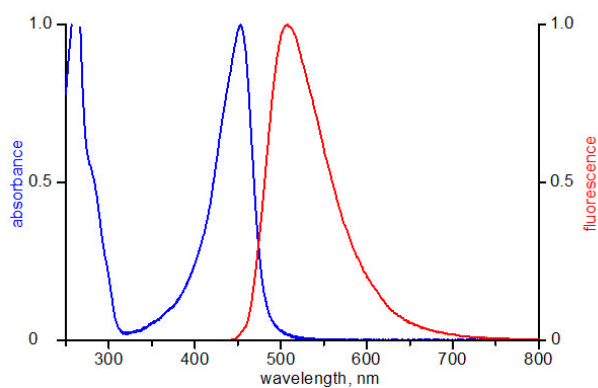
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 453 nm, λ_{em} 506 nm, ε 75.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ddATP-ATTO-465



excitation and emission spectrum of ATTO 465