



7-Propargylamino-7-deaza-ddATP-ATTO-390

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with ATTO 390, Triethylammonium salt

Cat. No.	Amount
NU-1612-390	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: C₃₄H₄₃N₆O₁₄P₃ (free acid)

Molecular Weight: 852.66 g/mol (free acid)

Exact Mass: 852.21 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

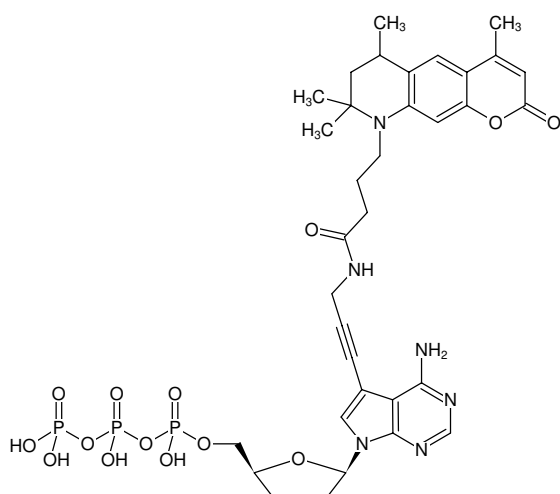
Form: solution in water

Color: slightly yellow

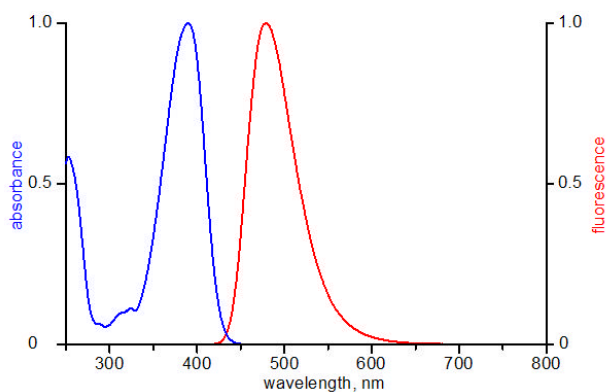
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 390 nm, λ_{em} 476 nm, ε 24.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ddATP-ATTO-390



excitation and emission spectrum of ATTO 390