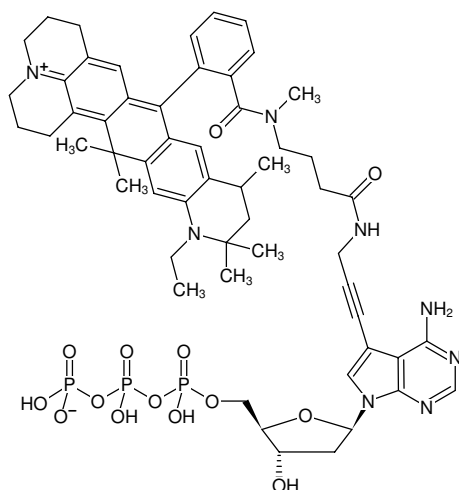




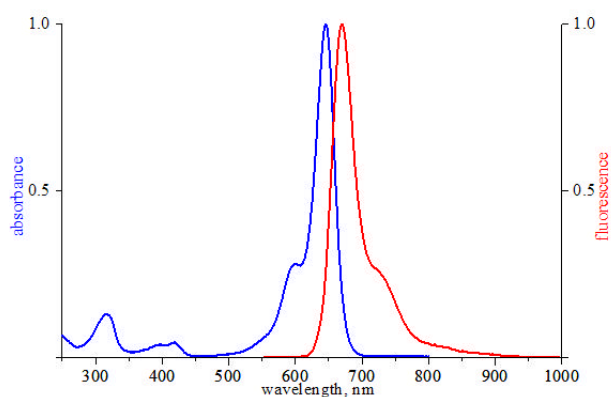
7-Propargylamino-7-deaza-dATP-ATTO-647N

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 647N, Triethylammonium salt

Cat. No.	Amount
NU-1611-647N	40 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dATP-ATTO-647N



excitation and emission spectrum of ATTO 647N

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: C₅₆H₆₉N₈O₁₄P₃ (free acid)

Molecular Weight: 1171.11 g/mol (free acid)

Exact Mass: 1170.41 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 646 nm, λ_{em} 664 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)