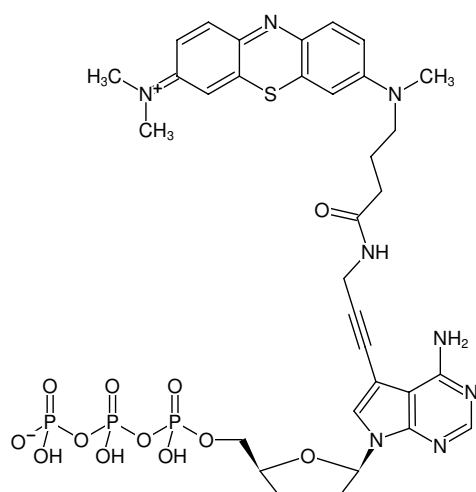




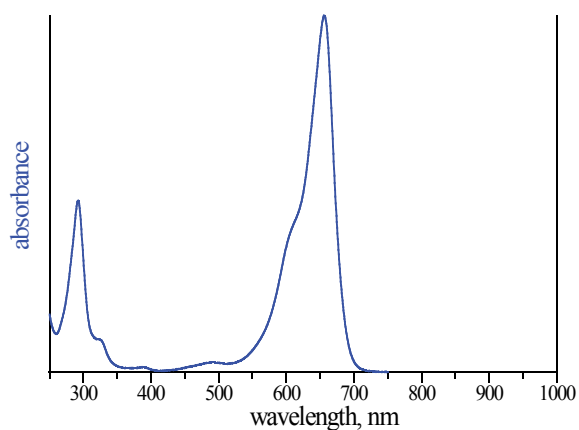
7-Propargylamino-7-deaza-ddATP-ATTO-MB2

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-1612-MB2	80 µl (0,5 mM)



Structural formula of 7-Propargylamino-7-deaza-ddATP-ATTO-MB2



absorption spectrum of ATTO MB2

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: C₃₃H₃₉N₈O₁₂P₃S (free acid)

Molecular Weight: 864.70 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{abs} 668 nm, ε 100.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)