



7-Propargylamino-7-deaza-ddGTP - ATTO-MB2

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

Cat. No.	Amount
NU-1618-MB2	80 µl (0,5 mM)

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months

Molecular Formula: C₃₃H₃₉N₈O₁₃P₃S (free acid)

Molecular Weight: 880.70 g/mol (free acid)

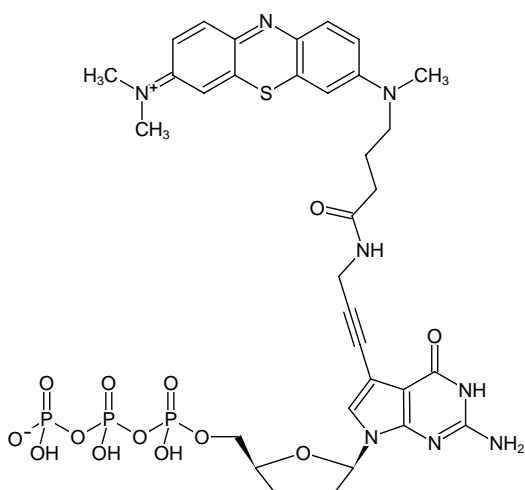
Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution, pH 7.5 ± 0.5

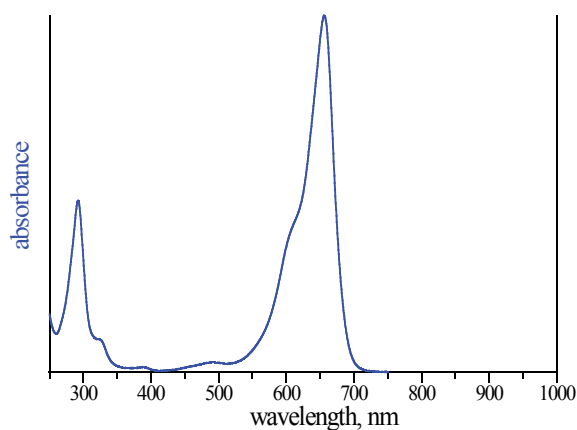
Concentration: 0.5 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{abs} 658 nm; ε 100.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



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



absorption spectrum of ATTO MB2