



5-Propargylamino-ddUTP-ATTO-MB2

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-1619-MB2	80 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: C₃₁H₃₇N₆O₁₄P₃S (free acid)

Molecular Weight: 842.64 g/mol (free acid)

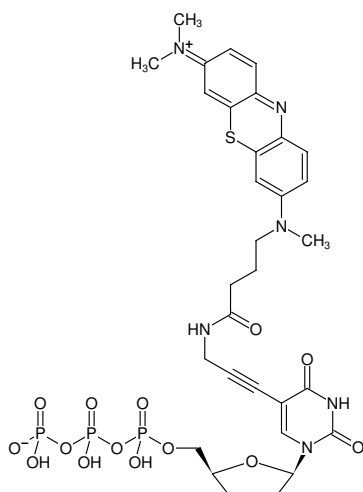
Purity: ≥ 95 % (HPLC)

Form: solution in water

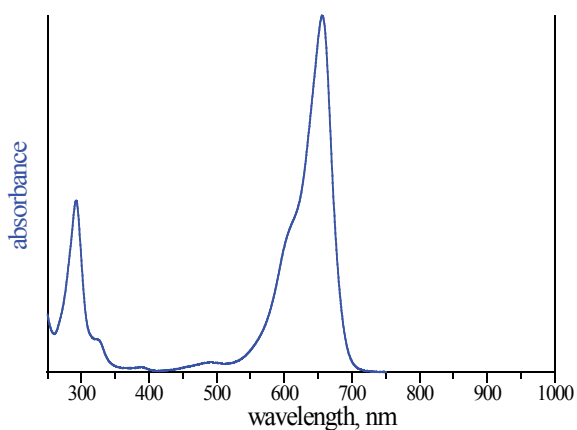
Concentration: 0.50 mM - 0.55 mM

pH: 7.5 ± 0.5

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



Structural formula of 5-Propargylamino-ddUTP-ATTO-MB2



absorption spectrum of ATTO MB2