

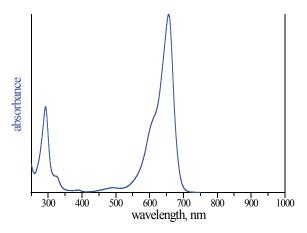


7-Propargylamino-7-deaza-ddGTP-ATTO-MB2

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)

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



absorption spectrum of ATTO MB2

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:** $C_{33}H_{39}N_8O_{13}P_3S$ (free acid) **Molecular Weight:** 880.70 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 $^{\text{-}1}$ cm $^{\text{-}1}$

(Tris-HCl pH 7.5)