



7-Propargylamino-7-deaza-ddATP-DYQ-661

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with DYQ 661, Triethylammonium salt

Cat. No.	Amount
NU-1612-661Q	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.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₃H₆₁N₇O₁₉P₃S₂ (Anion)

Molecular Weight: 1257.14 g/mol (Anion)

Purity: ≥ 95 % (HPLC)

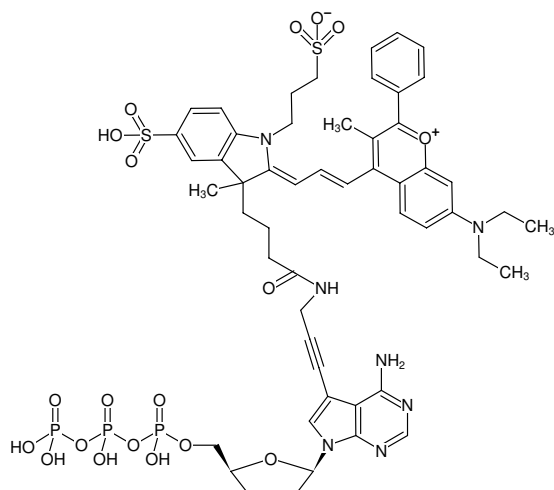
Form: clear aqueous solution

Concentration: 0.50 mM - 0.55 mM

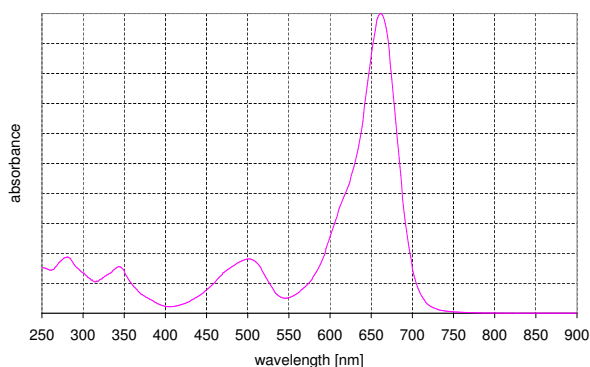
pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 662 nm, ε 104.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

λ_{exc} 662 nm, ε 140.0 L mmol⁻¹ cm⁻¹ (Ethanol)



Structural formula of 7-Propargylamino-7-deaza-ddATP-DYQ-661



absorption spectrum of DYQ 661

Please note: DYQ661 is not a dark quencher, λ_{em} 690 nm^[1]

Selected References:

[1] Tan *et al.* (2010) Direct detection of Delta9-tetrahydrocannabinol in saliva using a novel homogeneous competitive immunoassay with fluorescence quenching. *Anal. Chim. Acta.* **658** (2):187.