



7-Propargylamino-7-deaza-dGTP-DY-776

7-Deaza-7-propargylamino-2'-deoxyguanosine-5'-triphosphate, labeled with DY 776, Triethylammonium salt

Cat. No.	Amount
NU-1615-776	40 μ l (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume \leq 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: C₅₈H₆₆N₇O₂₁P₃S₂ (free acid)

Molecular Weight: 1354.23 g/mol (free acid)

Purity: \geq 95 % (HPLC)

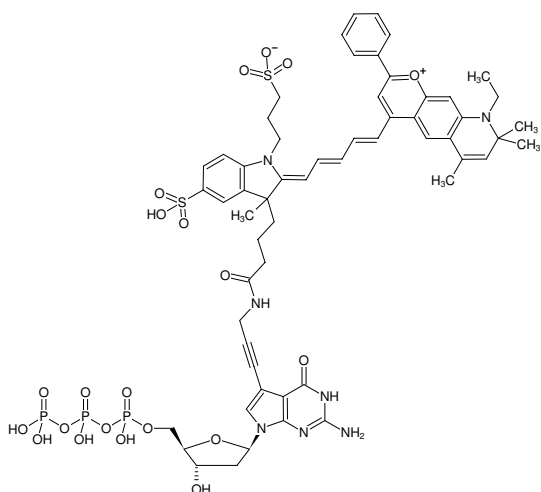
Form: solution in water

Color: dark blue

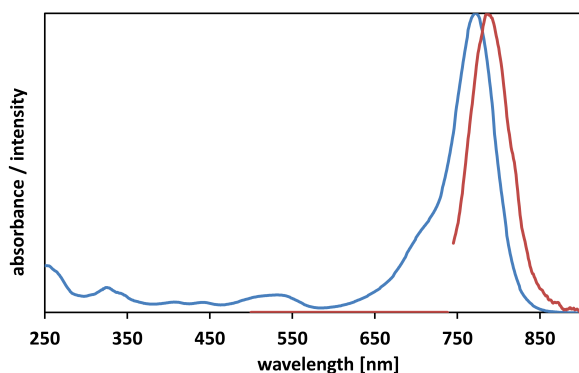
Concentration: 1.0 mM - 1.1 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{exc} 771 nm, λ_{em} 793 nm, ϵ 240.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-dGTP-DY-776



excitation and emission spectrum of DY 776