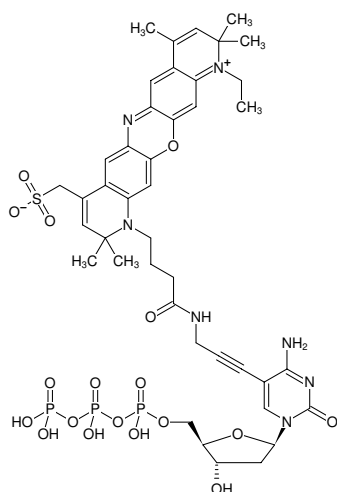




5-Propargylamino-dCTP-ATTO-700

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with ATTO 700, Triethylammonium salt

Cat. No.	Amount
NU-809-700	40 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-ATTO-700

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: C₄₂H₅₂N₇O₁₈P₃S (free acid)

Molecular Weight: 1067.89 g/mol (free acid)

Exact Mass: 1067.23 g/mol (free acid)

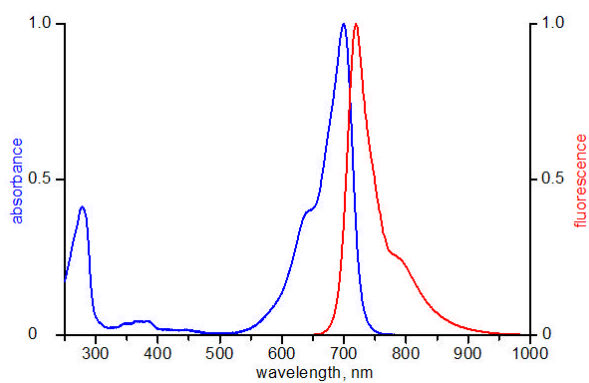
Purity: ≥ 95 % (HPLC)

Form: solution in water

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 700 nm, λ_{em} 716 nm, ε 120.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)



excitation and emission spectrum of ATTO 700