

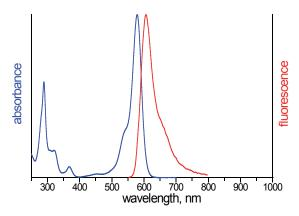


5-Propargylamino-dCTP-ATTO-Thio12

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with ATTO Thio12, Triethylammonium salt

Cat. No.	Amount
NU-809-THIO12	40 μl (1 mM)

Structural formula of 5-Propargylamino-dCTP-ATTO-Thio12



excitation and emission spectrum of ATTO Thio12

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 $\textbf{Molecular Formula:} \ C_{41}H_{48}N_7O_{15}P_3S \ (free \ acid)$ Molecular Weight: 1003.84 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} 582 nm, λ_{em} 607 nm, ϵ 110.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)