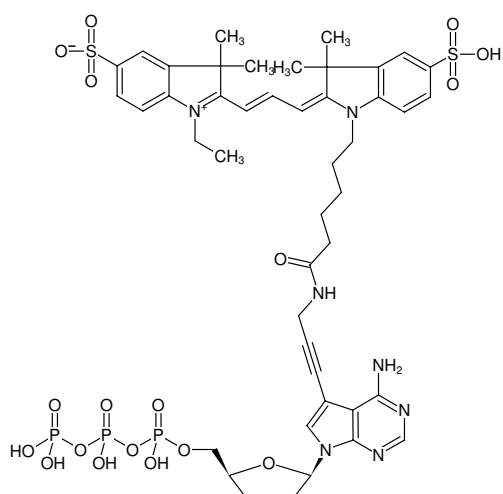




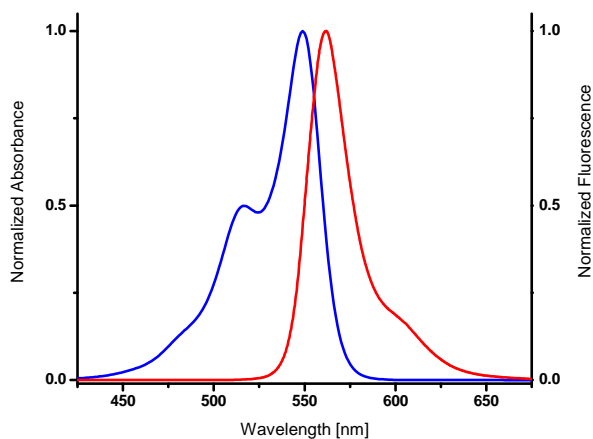
7-Propargylamino-7-deaza-ddATP-Cy3

7-Deaza-7-propargylamino-2',3'-dideoxyadenosine-5'-triphosphate, labeled with Cy3, Triethylammonium salt

Cat. No.	Amount
NU-1612-CY3	30 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-ddATP-Cy3



excitation and emission spectrum of Cy3

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: C₄₅H₅₆N₇O₁₈P₃S₂ (free acid)

Molecular Weight: 1140.01 g/mol (free acid)

Exact Mass: 1139.23 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: pink

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 550 nm, λ_{em} 570 nm, ε 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Masimba *et al.* (2014) Development of a simple microarray for genotyping HIV-1 drug resistance mutations in the reverse transcriptase gene in rural Tanzania. *Trop. Med. Int. Health.* **19** (6):664.

Cramer *et al.* (2007) Rapid microarray-based method for monitoring of all currently known single-nucleotide polymorphisms associated with parasite resistance to antimalaria drugs. *J. Clin. Microbiol.* **45** (11):3685.