



7-Propargylamino-7-deaza-dATP-Cy5

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-1611-CY5-S	25 µl (1 mM)
NU-1611-CY5-L	5 x 25 µl (1 mM)

For general laboratory use.

Please centrifuge briefly before opening (volume ≤ 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: C₄₇H₅₆N₇O₁₉P₃S₂ (free acid)

Molecular Weight: 1182.05 g/mol (free acid)

Exact Mass: 1181.24 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in 10 mM Tris-HCl

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{exc} 649 nm, λ_{em} 670 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

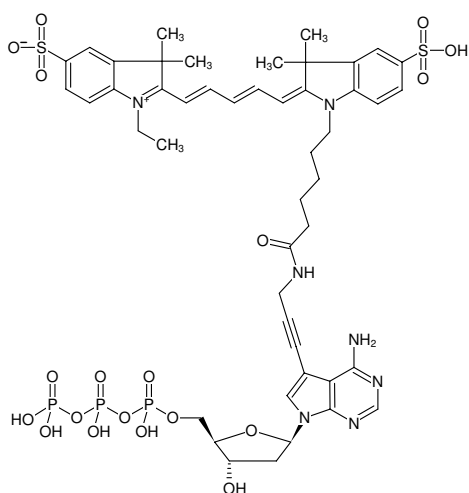
Incorporation by primer extension^[1]

Incorporation by PCR^[2]

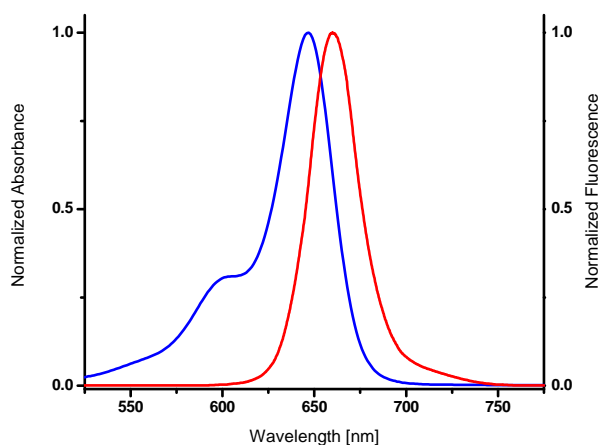
Selected References:

[1] Braslavsky *et al.* (2003) Sequence information can be obtained from single DNA molecules. *Proc. Natl. Acad. Sci. USA.* **100** (7):3960.

[2] Tasara *et al.* (2003) Incorporation of reporter molecule-labeled nucleotides by DNA polymerases. II. High-density labeling of natural DNA. *Nucleic Acids Res.* **31** (10):2636.



Structural formula of 7-Propargylamino-7-deaza-dATP-Cy5



excitation and emission spectrum of Cy5