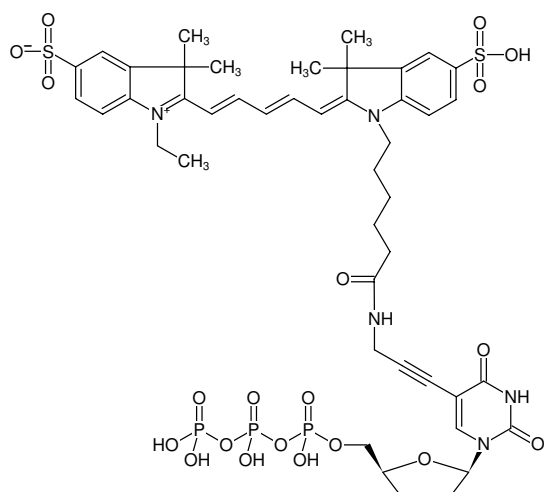




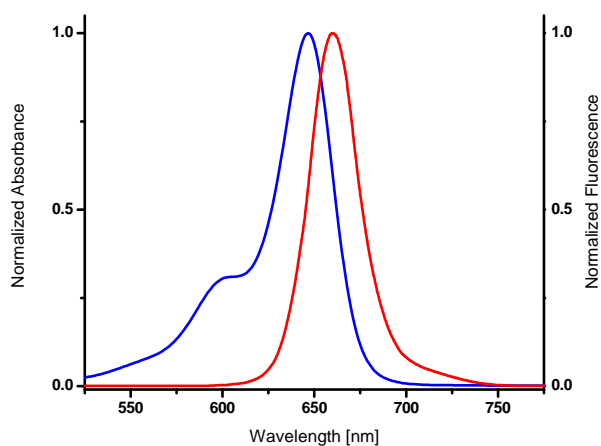
## 5-Propargylamino-ddUTP-Cy5

5-Propargylamino-2',3'-dideoxyuridine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-1619-CY5	30 µl (1 mM)



Structural formula of 5-Propargylamino-ddUTP-Cy5



excitation and emission spectrum of Cy5

**For research use only!**

**Shipping:** shipped on blue ice

**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:** C<sub>45</sub>H<sub>56</sub>N<sub>5</sub>O<sub>20</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1144.00 g/mol (free acid)

**Exact Mass:** 1143.22 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** blue

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 649 nm, λ<sub>em</sub> 670 nm, ε 250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.

Fredriksson *et al.* (2004) Assessing hematopoietic chimerism after allogeneic stem cell transplantation by multiplexed SNP genotyping using microarrays and quantitative analysis of SNP alleles. *Leukemia* **18** (2):255.

van Moorsel *et al.* (2004) beta-Globin mutation detection by tagged single-base extension and hybridization to universal glass and flow-through microarrays. *Eur. J. Hum. Genet.* **12** (7):567.