



### 5-Propargylamino-ddCTP-5/6-TAMRA

5-Propargylamino-2',3'-dideoxycytidine-5'-triphosphate, labeled with 5/6-TAMRA, Triethylammonium salt

Cat. No.	Amount
NU-850-TAM	30 µl (1 mM)

**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:** C<sub>37</sub>H<sub>39</sub>N<sub>6</sub>O<sub>16</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 916.66 g/mol (free acid)

**Exact Mass:** 916.16 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** pink to red

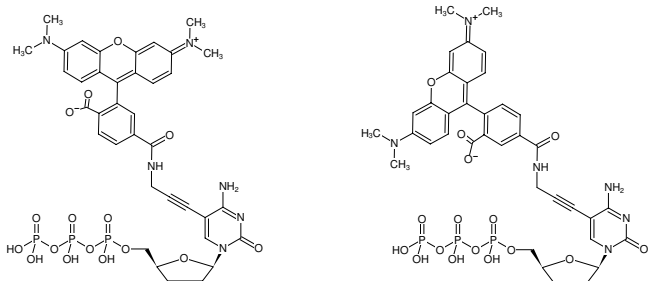
**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

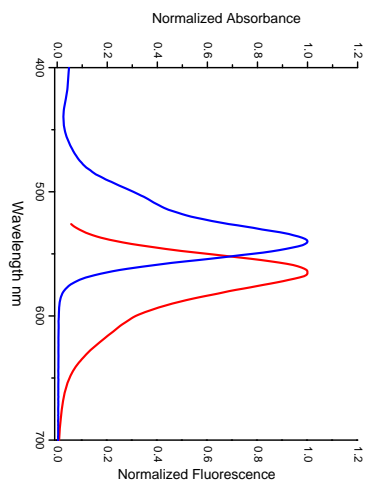
**Spectroscopic Properties:** λ<sub>exc</sub> 545 nm, λ<sub>em</sub> 575 nm, ε 90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

Esteves *et al.* (2011) Clinical relevance of multiple single-nucleotide polymorphisms in *Pneumocystis jirovecii* Pneumonia: development of a multiplex PCR-single-base-extension methodology. *J. Clin. Microbiol.* **49** (5):1810.



Structural formula of 5-Propargylamino-ddCTP-5/6-TAMRA



excitation and emission spectrum of 5/6-TAMRA