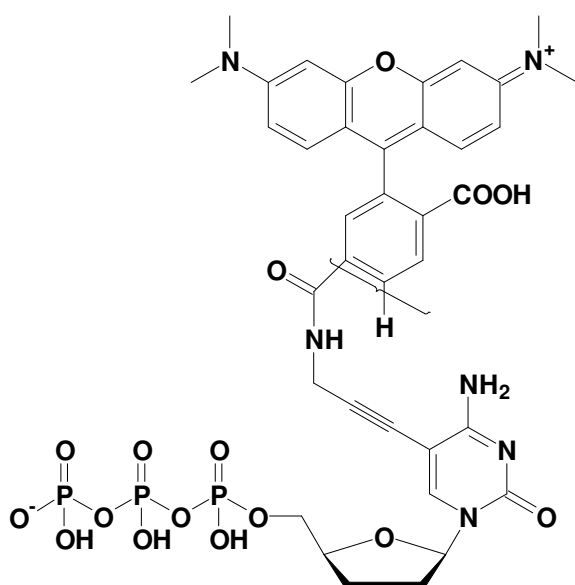


5-Propargylamino-ddCTP - 5/6-TAMRA

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

Cat. No.	Amount
NU-850-TAM	60 μ l / 0,5 mM



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

Cat. No.: NU-850-TAM

Molecular Formula: C₃₇H₃₉N₆O₁₆P₃ (free acid)

Molecular Weight: 916.66 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{exc} 545 nm; λ_{em} 575 nm; ϵ 90,000 cm⁻¹ M⁻¹

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

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.