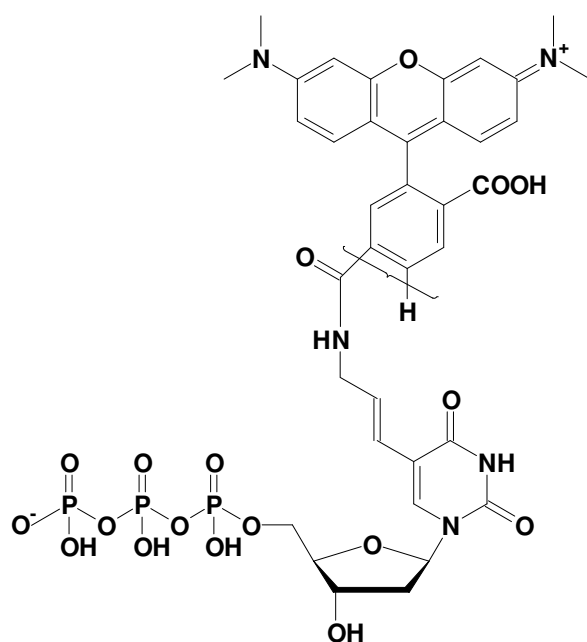


Aminoallyl-dUTP - 5/6-TAMRA

5-(3-Aminoallyl)-2'-deoxy-uridine-5'triphosphate labeled with 5/6-TAMRA, Triethylammonium salt

Cat. No.	Amount
NU-803-TAM	30 μ l (1 mM)



structural formula of Aminoallyl-dUTP - 5/6-TAMRA

Molecular Formula: $C_{37}H_{40}N_5O_{18}P_3$ (free acid)

Molecular Weight: 935.66 g/mol (free acid)

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

Spectroscopic properties: λ_{abs} 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!