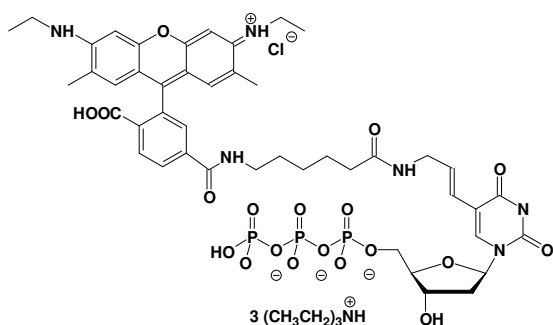


Rhodamin-12-dUTP

Rhodamin-X-5-Aminoallyl-2'-desoxy-uridine-5'-triphosphate, triethylammonium salt

Cat. No.	Amount
NU-803-RHOX	30 μ l / 1 mM



Cat. No.: NU-803-RHOX

Molecular Formula: C₃₉H₄₃N₆O₁₉P₃ (Anion)

Molecular Weight: 990.70 (Anion)

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

Spectroscopic properties:

λ_{exc} 505 nm; λ_{em} 530 nm; ϵ 85,000 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!