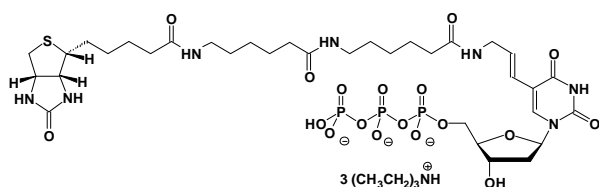


Biotin-18-dUTP

γ -[N-(Biotin-6-amino-hexanoyl-6-aminohexanoyl)]-5-(3-aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-BIOXX	10 μ l / 5 mM



Cat. No.: NU-803-BIOXX

Molecular Formula: C₃₄H₅₃N₇O₁₈P₃S (Anion)

Molecular Weight: 972.81 (Anion)

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

Spectroscopic properties: λ_{\max} 240 nm; ϵ 10700

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!