

Biotin-18-dCTP

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

Cat. No.	Amount
NU-809-BIOXX	200 μ l / 5 mM

Cat. No.: NU-809-BIOXX

Molecular Formula: C₃₄H₅₅N₈O₁₇P₃S (free acid)

Molecular Weight: 972.83 (free acid)

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

Spectroscopic properties: λ_{\max} 294 nm;
 ϵ 9300

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!