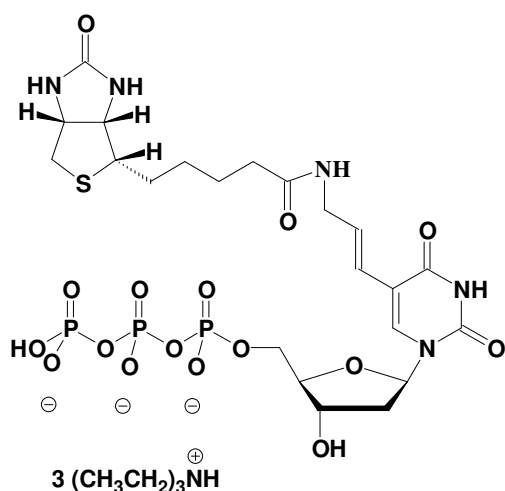


Biotin-dUTP

N-Biotin-5-(3-aminoallyl)-2'-deoxy-uridine-5'triphosphate, Triethylammonium salt

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



Cat. No.: NU-803-BIO

Molecular Formula: C₂₂H₃₃N₅O₁₆P₃S (Anion)

Molecular Weight: 748.52 (Anion)

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

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!