

Biotin-4-UTP

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

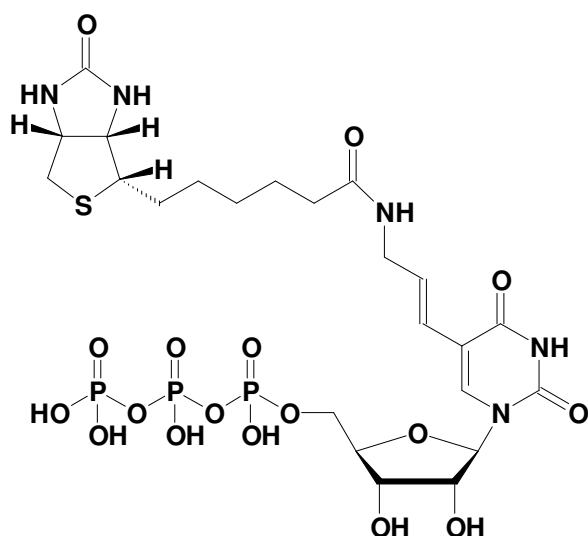
Cat. No.	Amount
NU-821-BIO	60 μ l / 5 mM

Applications:

RNA-labeling^[1]

Selected References:

[1] Fen *et al.* (1990) Direct quantitation of biotin-labeled nucleotide analogs in RNA transcripts. *Analytic. Biochem.* **190**:78.



Cat. No.: NU-821-BIO

Molecular Formula: C₂₃H₃₆N₅O₁₇P₃S (free acid)

Molecular Weight: 779.54 (free acid)

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

Spectroscopic properties: λ_{\max} 249/289 nm;
 ϵ 10700/7100

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!