

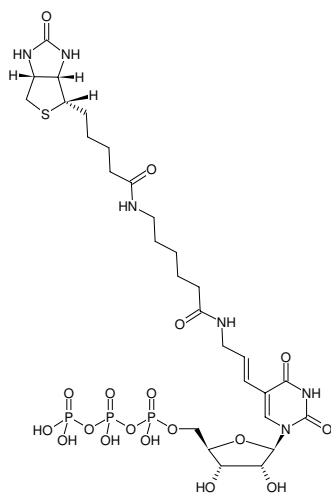


Biotin-11-UTP - high concentration

Biotin-X-5-Aminoallyl-UTP

Biotin-X-5-Aminoallyl-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-BIOX-HC	30 µl (75 mM)



Structural formula of Biotin-11-UTP - high concentration

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₂₈H₄₅N₆O₁₈P₃S (free acid)

Molecular Weight: 878.67 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: sterile clear aqueous solution in 20 mM Tris-HCl

Concentration: 75 mM - 80 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 240/289 nm, ε 10.7/71 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Microarray based hybridization detection^[1, 2]

RNA-labeling^[2]

Biotin-PAGE^[3]

ISH^[4]

Selected References:

[1] Lonergan *et al.* (2007) Comparison of target labeling methods for use of Affymetrix GeneChips. *BMC Biotechnology* **7**:24.

[2] Glaid *et al.* (1989) Non-isotopic RNA probes. Comparison between different labels and detection systems. *Histochemistry* **93**:91.

[3] Theissen *et al.* (1989) Degree of biotinylation in nucleic acids estimated by a gel retardation assay. *Analyt. Biochem.* **179**:98.

[4] Baumann *et al.* (1988) Flow cytometric detection of ribosomal RNA in suspended cells by fluorescent in situ hybridization. *Cytometry* **9**:517.