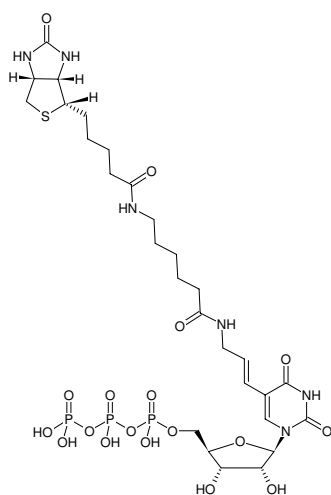


**Biotin-11-UTP**

Biotin-X-(5-aminoallyl)-UTP

Biotin-X-(5-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-BIOX	30 µl (10 mM)



Structural formula of Biotin-11-UTP

For research use only!**Shipping:** shipped on gel packs**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)**Exact Mass:** 878.17 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** filtered solution in 10 mM Tris-HCl**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 289 nm, ε 7.1 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]**Related Products:**

HighYield T7 RNA Synthesis Kit, #RNT-101

HighYield T7 Biotin11 RNA Labeling Kit, #RNT-101-BIOX

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.