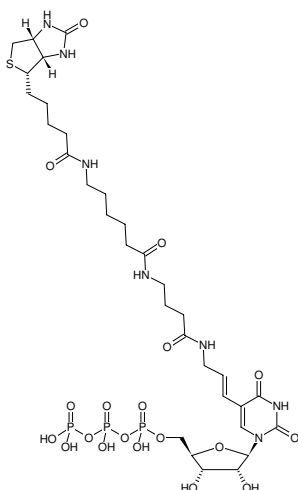


**Biotin-16-UTP**

Biotin-16-(5-aminoallyl)-UTP

Biotinyl-ε-aminocaproyl-γ-aminobutyryl-5-(3-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-BIO16	30 µl (10 mM)



Structural formula of Biotin-16-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:** C₃₂H₅₂N₇O₁₉P₃S (free acid)**Molecular Weight:** 963.78 g/mol (free acid)**Exact Mass:** 963.23 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} 240/289 nm, ε 10.7/7.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**Microarray based hybridization detection^[1,2]Northern Blot^[3]**Related Products:**

HighYield T7 RNA Synthesis Kit, #RNT-101

HighYield T7 Biotin16 RNA Labeling Kit, #RNT-101-BIO16

Selected References:

[1] Li et al. (2006) Identification of estrogen-responsive genes in parenchyma and fat pad of the bovine mammary gland by microarray analysis. *Physiol. Genomics* **27**:42.

[2] Kim et al. (2007) Galanin is up-regulated in colon adenocarcinoma. *Cancer Epidemiol. Biomarkers Prev.* **16**:2373.

[3] Jelinsky et al. (1999) Global response of *Saccharomyces cerevisiae* to alkylating agent. *Proc. Natl. Acad. Sci. USA* **96**:1486.