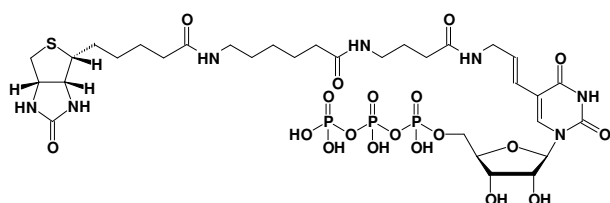


## Biotin-16-UTP

$\gamma$ -[N-(Biotin-6-amino-hexanoyl-6-aminobutanoyl)]-5-(3-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-821-BIO16	60 $\mu$ l / 5 mM



**Cat. No.:** NU-821-BIO16

**Molecular Formula:** C<sub>32</sub>H<sub>52</sub>N<sub>7</sub>O<sub>19</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 963.78 (free acid)

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

**Spectroscopic properties:**  $\lambda_{\max}$  240 nm;  $\epsilon$  10,700 M<sup>-1</sup>cm<sup>-1</sup>

### 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!**

### Applications:

Microarray based hybridization detection<sup>[1, 2]</sup>

Northern Blot<sup>[3]</sup>

### 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.