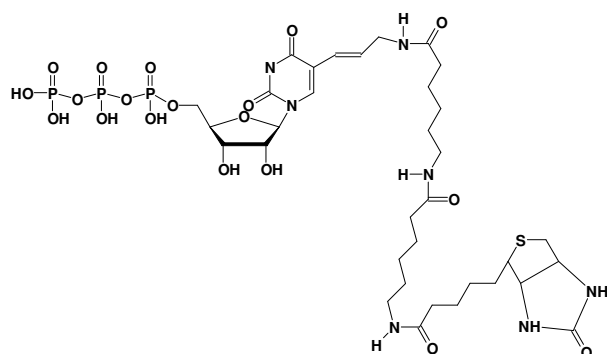


Biotin-18-UTP

γ -[N-(Biotin-6-amino-hexanoyl-6-aminohexanoyl)]-5-(3-aminoallyl)-uridine-5'-triphosphate, Triethylammonium salt

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



allowing simultaneous visualization by the use of confocal microscopy of three cellular mRNA species in individual neurons. *J Histochem Cytochem.* **46(6)**:753.

Strissel *et al.* (1996) Scaffold attachment regions in centromere-associated DNA. *Chromosoma.* **105(2)**:122.

Cat. No.: NU-821-BIOXX

Molecular Formula: C₃₃H₅₃N₇O₁₉P₃S (free acid)

Molecular Weight: 976.80 (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!

Selected References:

Stoffels *et al.* (1999) rRNA probe-based cell fishing of bacteria. *Environ Microbiol.* **1(3)**:259.

Grino *et al.* (1998) An in situ hybridization histochemistry technique