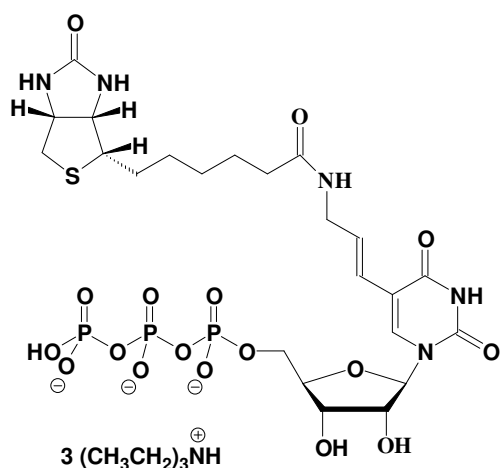


## Biotin-UTP

$\gamma$ -N-(Biotin)-5-(3-aminoallyl)-uridine-5'triphosphate, Triethylammonium salt

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



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

**Molecular Formula:**  $C_{23}H_{33}N_5O_{17}P_3S$  (Anion)

**Molecular Weight:** 776.51 (Anion)

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

**Spectroscopic properties:**  $\lambda_{max}$  249/289 nm;  
 $\epsilon$  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!**