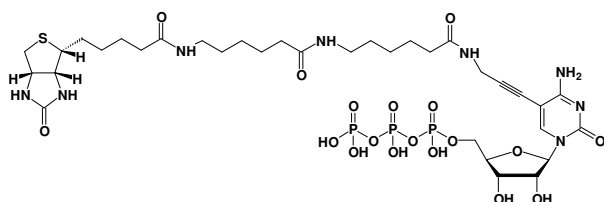


## Biotin-18-CTP

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

| Cat. No.     | Amount            |
|--------------|-------------------|
| NU-831-BIOXX | 60 $\mu$ l / 5 mM |



**Cat. No.:** NU-831-BIOXX

**Molecular Formula:**  $C_{34}H_{55}N_8O_{18}P_3S$  (free acid)

**Molecular Weight:** 988.83 (free acid)

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

**Spectroscopic properties:**  $\lambda_{max}$  294 nm;  
 $\epsilon$  9300

### 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!**