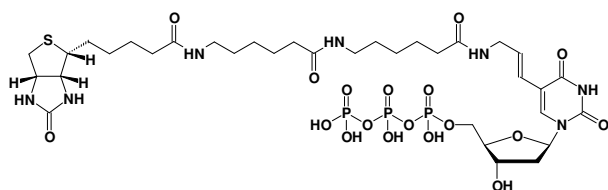


## Biotin-18-dUTP

$\gamma$ -[N-(Biotin-6-amino-hexanoyl-6-aminohexanoyl)]-5-(3-aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-BIOXX	200 $\mu$ l / 5 mM



**Cat. No.:** NU-803-BIOXX

**Molecular Formula:** C<sub>34</sub>H<sub>56</sub>N<sub>7</sub>O<sub>18</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 975.83 (free acid)

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

**Spectroscopic properties:**  $\lambda_{\max}$  240 nm;  $\epsilon$  10700

### 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!**