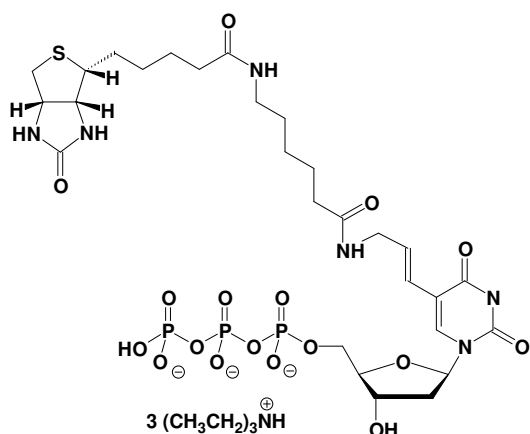


Biotin-11-dUTP

Biotin-X-5-Aminoallyl-2'-desoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-BIOX	10 μ l / 5 mM



Cat. No.: NU-803-BIOX

Molecular Formula: $C_{28}H_{41}N_6O_{17}P_3S$ (Anion)

Molecular Weight: 858.64 (Anion)

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

Spectroscopic properties: λ_{max} 240/289 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!