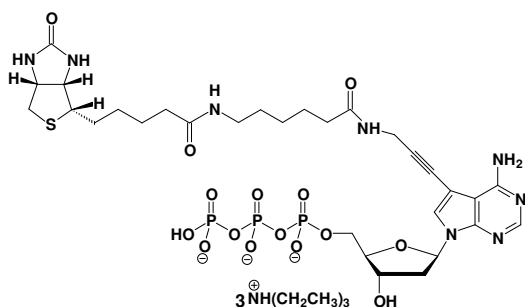


Biotin-11-dATP

γ -[N-(Biotin-6-amino-hexanoyl)]-7-propargylamino-2'-deoxy-7-deaza-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1175-BIOX	10 μ l / 5 mM



Cat. No.: NU-1175-BIO

Molecular Formula: $C_{30}H_{42}N_8O_{15}P_3S$ (Anion)

Molecular Weight: 879.68 (Anion)

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

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!