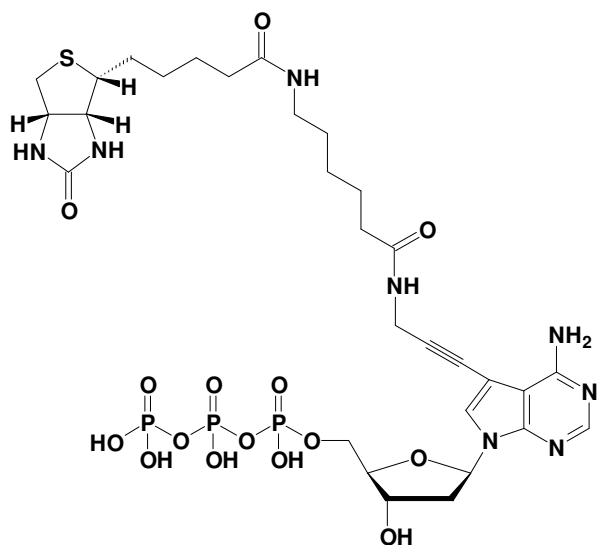


Biotin-11-dATP

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: C₃₀H₄₅N₈O₁₅P₃S (free acid)

Molecular Weight: 882.71 (free acid)

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

Spectroscopic properties: λ_{\max} 280 nm;
 ϵ 12,700

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!

Selected References:

Vorwerk *et al.* (2008) Microfluidic-based enzymatic on-chip labeling of miRNAs. *N. Biotechnol.* **25** (2-3):142.

Zammatteo *et al.* (2005) Unambiguous identification of the expressed MAGE-A genes on a DNA microarray. *Clin. Chem.* **51** (12):2420.

Dauwerse *et al.* (1999) Two-colour FISH detection of the inv(16) in interphase nuclei of patients with acute myeloid leukaemia. *Br. J. Haematol.* **106** (1):111.

Wright *et al.* (1998) Proliferating and migrating mesangial cells responding to injury express a novel receptor protein-tyrosine phosphatase in experimental mesangial proliferative glomerulonephritis. *J. Biol. Chem.* **273** (37):23929.