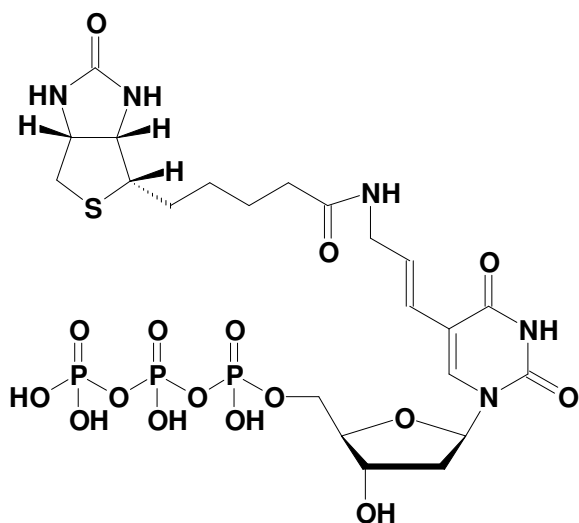


Biotin-4-dUTP

N-Biotin-5-(3-aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-BIO	200 μ l / 5 mM



Cat. No.: NU-803-BIO

Molecular Formula: C₂₂H₃₄N₅O₁₆P₃S (free acid)

Molecular Weight: 749.51 (free acid)

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!

Applications:

In situ hybridization^[1]

SNP-Analysis^[2, 3]

Excision repair patches^[4]

Southern Blot^[5, 6]

DNA-labeling^[7, 8]

Selected References:

[1] Baechner *et al.* (1999) Apolipoprotein E, a Bmp-2 upregulated gene in mesenchymal progenitors, is highly expressed in murine embryonic development. *Biofactors* **9**:11.

[2] Yanagawa *et al.* (2009) Modified multiple primer extension method. *Methods in Molecular Biology* **578**:425.

[3] Litos *et al.* (2009) Multianalyte, dipstick-type, nanoparticle-based DNA biosensors for visual genotyping of single-nucleotide polymorphisms. *Biosensors and Bioelectronics* **24**:3135.

[4] Larone *et al.* (1991) Purification of DNA fragments containing excision-repair patches from human cells using streptavidin-biotin. *Mutation Research, DNA Repair* **254**:273.

[5] Lauc *et al.* (1994) A non-radioactive, sensitive method for detection of DNA fragmentation in apoptotic cells. *Anal. Cell Pathol.* **2**:107.

[6] Gregersen *et al.* (1987) Improved methods for the detection of unique sequences in Southern blots of mammalian DNA by non-radioactive biotinylated DNA hybridization probes. *Clinica Chimica Acta* **169**:267.

[7] Langer *et al.* (1981) Enzymatic synthesis of biotin-labeled polynucleotides: Novel nucleic acid affinity probes. *Proc. Natl. Acad. Sci. USA* **78**:6633.

[8] Tasara *et al.* (2003) Incorporation of reporter molecule-labeled nucleotides by DNA polymerases. II. High density labeling of natural DNA. *Nucleic Acids Res.* **31(10)**:2636.