



Biotin-11-dUTP, 1 mM in TE buffer

Catalog Number	Packaging Size
C403A	50 µL

Storage upon receipt: -20°C

User Manual

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, MD 20850

USA

Tel: 1-866-360-9531; 1-301-762-0888

Fax: 1-301-762-3888

Inquiry: inquiry@genecopoeia.com

Order: orders@genecopoeia.com

Web: www.genecopoeia.com

I. Product Description

Biotin-11-dUTP (biotin- ϵ -aminocaproyl-[5-{3-aminoallyl}-2'-deoxyuridine-5'-triphosphate]) is supplied as 1 mM solution in TE buffer. The nucleotide is designed for enzymatic non-radioactive labeling of DNA by PCR, nick-translation, cDNA synthesis, random primed labeling, or primer extension. Biotin-11-dUTP can be enzymatically incorporated into DNA with Reverse Transcriptases, Taq DNA Polymerase, phi29 DNA Polymerase, Klenow Fragment, exo-, Klenow Fragment, and DNA Polymerase I. The number '11' is the number of atoms in the linker between biotin and dUTP.

II. Specifications

Concentration: 1 mM solution in TE buffer

Storage: Store at -20°C

Molecular Formula: $C_{28}H_{42}N_6Na_3O_{17}P_3S$

Molecular Weight: 928.6

III. General Characteristics

$\lambda_{max}=240\text{ nm}$, $\epsilon=10.7 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$ (pH 7.0);

$\lambda_{max}=289\text{ nm}$, $\epsilon = 7.1 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$ (pH 7.0).

