



## Biotin-11-UTP, 1 mM in TE buffer

Catalog Number	Packaging Size
C404A	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](mailto:inquiry@genecopoeia.com)

Order: [orders@genecopoeia.com](mailto:orders@genecopoeia.com)

Web: [www.genecopoeia.com](http://www.genecopoeia.com)

## I. Product Description

Biotin-11-UTP is supplied as 1 mM solution in TE buffer. The nucleotide is designed for indirect nonradioactive enzymatic labeling of RNA during *in vitro* transcription. Biotin-11-UTP can be enzymatically incorporated into RNA with T7, T3 and SP6 RNA polymerases, as well as other RNA modifying enzymes. The number '11' is the number of atoms in the linker between biotin and UTP. Biotinylated RNA can be used in place of radioactively labeled RNA in many applications, including cRNA amplification and labeling for microarray target preparation. Detection is then carried out via one of a variety of streptavidin-based methods.

## II. Specifications

Concentration: 1 mM solution in TE buffer

Storage: Store at -20°C

Molecular Formula:  $C_{28}H_{42}N_6Na_3O_{18}P_3S$

Molecular Weight: 944.62

## 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).

