



5-Aminoallyl-UTP (AA-UTP), lyophilized powder

Catalog Number	Packaging Size
C402B	1 μ mol

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

5-Aminoallyl-UTP (5-[3-aminoallyl]-2'-uridine-5'-triphosphate) is supplied as lyophilized form. The nucleotide is designed for indirect nonradioactive enzymatic labeling of RNA during *in vitro* transcription. 5-Aminoallyl-UTP can be enzymatically incorporated into RNA with T7, T3 and SP6 RNA polymerases. The resulting amine-containing RNA can be subsequently labeled with any amine-reactive fluorescent dye, biotin, or hapten. This two-step method for labeling nucleic acids is considerably more economical than the one-step method using a prelabeled UTP.

II. Specifications

Concentration: NA

Storage: Store at -20°C

Molecular Formula: $C_{12}H_{17}N_3Na_3O_{15}P_3$

Molecular Weight: 605.17

III. General Characteristics

$\lambda_{max}=240\text{ nm}$, $\epsilon=11.9 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$ (pH 7.0);

$\lambda_{max}=290\text{ nm}$, $\epsilon=7.8 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$ (pH 7.0).

