



## Cy3-X-UTP, lyophilized powder

Catalog Number	Packaging Size
C420B	25 nmol

**Storage upon receipt:** -20°C, protected from light.

### 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

Cy3-X-UTP is supplied as lyophilized powder. The nucleotide is designed for indirect nonradioactive enzymatic labeling of RNA during *in vitro* transcription. Cy3-X-UTP can be enzymatically incorporated into RNA with T7, T3 and SP6 RNA polymerases, as well as other RNA modifying enzymes. The X linker between Cy3 and UTP can improve the incorporation efficiency.

## II. Specifications

Excitation/Emission (nm): 550/565 nm

Concentration: 1 mM in TE buffer

Storage: Store at -20°C

## III. General Characteristics

$\lambda_{max}=550\text{ nm}$ ,  $\epsilon=150.0 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$ .