



Andy Fluor 568-X-UTP, 1 mM in TE buffer

Catalog Number	Packaging Size
C414A	25 μ L

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

Order: orders@genecopoeia.com

Web: www.genecopoeia.com

I. Product Description

Andy Fluor 568-X-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. Andy Fluor 568-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 Andy Fluor 568 and UTP can improve the incorporation efficiency.

II. Specifications

Excitation/Emission (nm): 578/602 nm

Concentration: 1 mM in TE buffer

Storage: Store at -20°C

III. General Characteristics

$\lambda_{max}=578 \text{ nm}$, $\epsilon=91.0 \times 10^3 \text{ M}^{-1}\text{cm}^{-1}$.