



Andy Fluor 647-X-dUTP, 1 mM in TE buffer

Catalog Number	Packaging Size
C417A	25 µL

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 647-X-dUTP 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. Andy Fluor 647-X-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 X linker between Andy Fluor 647 and dUTP can improve the incorporation efficiency.

II. Specifications

Excitation/Emission (nm): 650/665 nm

Concentration: 1 mM in TE buffer

Storage: Store at -20°C

III. General Characteristics

$\lambda_{max}=650\text{ nm}$, $\epsilon=250.0 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$.